

User Guide

# SMC® Human IFN-λ1 High Sensitivity Immunoassay Kit

## Microparticle Assay

Human IFN-λ1 Immunoassay Kit for the Quantitative Determination of IFN-λ1 in Human Serum or Plasma

**03-0213-00**

|                                                   |    |                                                 |    |
|---------------------------------------------------|----|-------------------------------------------------|----|
| Introduction .....                                | 2  | Assay Reading .....                             | 12 |
| Supplies .....                                    | 3  | To read on the SMCxPRO® Immunoassay System..... | 12 |
| Reagents Included with the Kit .....              | 3  | SMC® Assay Overview .....                       | 13 |
| Kit Storage.....                                  | 3  | Assay Characteristics.....                      | 14 |
| Additional Supplies Required (Not provided) ..... | 4  | Sensitivity .....                               | 14 |
| Assay Best Practices .....                        | 5  | Precision .....                                 | 14 |
| Precautions .....                                 | 6  | Cross-Reactivity .....                          | 14 |
| Hazard Labels .....                               | 6  | Spike Recovery .....                            | 15 |
| Assay Preparation .....                           | 8  | Graph of Typical Reference Curve ..             | 15 |
| Sample Preparation .....                          | 8  | Troubleshooting .....                           | 16 |
| Sample Dilution.....                              | 8  | Well Plate .....                                | 18 |
| Reagent Preparation .....                         | 8  | Terms of Sale .....                             | 19 |
| Initial Standard Stock Preparation .....          | 9  | Notice .....                                    | 20 |
| Standard Curve.....                               | 9  | Technical Assistance .....                      | 20 |
| Assay Procedure .....                             | 10 | Terms and Conditions of Sale .....              | 20 |
| Target Capture.....                               | 10 | Safety Data Sheets (SDS) .....                  | 20 |
| Post-Capture Wash .....                           | 10 | Contact Information.....                        | 20 |
| Detection Antibody Incubation....                 | 11 |                                                 |    |
| Post-Detection Wash.....                          | 11 |                                                 |    |
| Post-Detection Shake.....                         | 11 |                                                 |    |
| Final Aspiration .....                            | 11 |                                                 |    |
| Elution.....                                      | 12 |                                                 |    |

---

## Introduction

Interferon lambda 1 (IFN- $\lambda$ 1), also known as IL-29, is a type of cytokine belonging to the type III interferon family, which also includes IL-28A and IL-28B. These cytokines are recognized for their role in the immune response against viral infections and are structurally like members of the IL-10 family. IL-29 is primarily produced by dendritic cells and is involved in the modulation of antiviral activity, acting similarly to type I interferons, though it interacts with a distinct receptor complex. This cytokine is crucial for its antiviral properties, immune regulation, and potential therapeutic implications in managing viral infections and certain autoimmune disorders. The specific signaling through its receptor can activate various pathways leading to the induction of antiviral state and regulation of immune responses.

The SMC<sup>®</sup> Human IFN- $\lambda$ 1 High Sensitivity Immunoassay uses a quantitative fluorescent sandwich immunoassay technique to measure IFN- $\lambda$ 1 in Human serum or plasma samples. A capture antibody specific for Human IFN- $\lambda$ 1 has been pre-coated onto paramagnetic microparticles (beads). The user pipettes beads, standards, and samples into uncoated microplate wells. During incubation, the IFN- $\lambda$ 1 present in the sample binds to the capture antibody on the coated beads. Unbound molecules are washed away during the subsequent wash steps. Fluor-labeled detection antibody is added to each well and incubated. This detection antibody recognizes and binds to IFN- $\lambda$ 1 that has been captured onto the beads, thus completing the sandwich. Elution buffer is added to dissociate the protein sandwich, releasing the fluor-labeled antibodies. The eluted antibodies are transferred to a SMC<sup>®</sup> 384-well Read Plate. The plate is loaded into the SMCxPRO<sup>®</sup> System where the labeled molecules are detected and counted. The number of fluor-labeled detection antibodies counted is directly proportional to the amount of IFN- $\lambda$ 1 present in the sample. The amount of IFN- $\lambda$ 1 in unknown samples is interpolated from a standard curve.

## Supplies

The SMC® Human IFN- $\lambda$ 1 High Sensitivity Immunoassay Kit includes all reagents listed below; these components are lot matched and not intended to be used separately. Additional reagents and supplies are required to run this immunoassay, as listed in the next section, Additional Supplies Required (not provided).

This kit and all reagents supplied are for research use only.

### Reagents Included with the Kit

All items are shipped with a cold pack unless otherwise stated.

| Description                         | Storage Conditions | Packaging Details  | Component Number |
|-------------------------------------|--------------------|--------------------|------------------|
| Assay Buffer                        | 2-8 °C             | 2 x 20 mL          | 02-9950-00       |
| IFN- $\lambda$ 1 Coated Beads       | 2-8 °C             | 1 x 550 $\mu$ L    | 02-2213-00       |
| Standard Diluent                    | 2-8 °C             | 2 x 20 mL          | 02-0225-02       |
| IFN- $\lambda$ 1 Detection Antibody | 2-8 °C             | 1 x 270 $\mu$ L    | 02-1213-00       |
| IFN- $\lambda$ 1 Standard           | 2-8 °C             | 1 lyophilized vial | 02-8213-00       |
| 10X Wash Buffer                     | 2-8 °C             | 3 x 50 mL          | 02-0001-03       |
| Buffer D                            | 2-8 °C             | 1 x 6 mL           | 02-0446-00       |
| Elution Buffer B                    | 2-8 °C             | 1 x 5 mL           | 02-0211-02       |

### Kit Storage

The SMC® Human IFN- $\lambda$ 1 High Sensitivity Immunoassay Kit should be stored at 2-8 °C. Discard standards after one use.

Supplied 10X Wash Buffer does not contain preservative. After dilution, the 1X Wash Buffer may be filter sterilized with Stericup® filter, for storage of up to 1 month at 2-8 °C. If not filter sterilized, all remaining 1X wash buffer should be discarded upon experiment completion.

Proper kit performance can only be guaranteed if the materials are stored properly.

## Additional Supplies Required (Not provided)

Catalogue numbers are provided to purchase products at [SigmaAldrich.com](https://www.sigmaaldrich.com) or through sales quote, unless otherwise noted.

### Equipment

- SMCxPRO® Ultrasensitive Immunoassay System for sample acquisition (95-0100-00)
- Orbital microplate shaker for assay plate incubation (for example, Boekel Scientific Jitterbug™ Shaker)
- BioTek® 405 TSUVS Microplate Washer for assay plate washing (95-0004-05)
- Sphere Mag Plate for performing microparticle capture (90-0003-02)
- Rotisserie tube rotator for microparticle suspension
- Benchtop centrifuge with bucket rotors capable of reaching 1,100 x g for sample/plate centrifugation
- Microcentrifuge capable of reaching 13,000 x g for reagent/sample centrifugation
- Single channel manual pipettes to accurately dispense 10-20 µL and 20-250 µL
- 12-channel manual pipettes to accurately dispense 10-20 µL and 20-250 µL
- Plate roller for complete plate sealing (Fisher Scientific, NC9185793)

### Supplies

- Micro-centrifuge tubes for sample preparation and storage
- 2 L Container with cap for Wash Buffer dilution
- Stericup® Quick Release Vacuum Filtration System, 0.22 µm, 1 L; for filter sterilizing 1X Wash Buffer (S2GPU11RE)
- MultiScreen®<sub>HTS</sub> 96-well Plate, hydrophilic PVDF membrane (MSBVN1210)
- 15 mL conical tube with cap for capture bead and detection antibody dilution
- 96-well V-bottom plate for assay setup (AXYP96450VCS)
- Axygen™ Microplate Sealing Film and Tapes (Fisher Scientific, 14-222-344)
- Universal plate cover to minimize plate well contamination (Fisher Scientific, 253623)
- 12-Channel reagent reservoir (sterile) for standard serial dilution (Argos/Cole Parmer, 04395-33)
- VistaLab® 25 mL Reservoirs for addition of reagents (Fisher Scientific, 21-381-27C)
- Millex® Syringe Filter, 0.2 µm for detection antibody filtration (SLGPR33RS)
- Luer-Lok® Syringe, 5 mL; for Detection Antibody Filtration (Fisher Scientific, 14-829-45)
- SMC® 384-well plate with adhesive seal (02-1008-00)
- SMC® 384-well plate, bulk case of 32 (ABB2-00160A)
- Nunc™ Aluminum adhesive plate seals (Fisher Scientific, 276014)

### Reagents

- 10X Wash Buffer for automated assay plate washing, 1 L (02-0111-00)
- De-ionized or distilled water for dilution of 10X Wash Buffer

For research use only. Not for use in diagnostic procedures.

## Assay Best Practices

To obtain reliable and reproducible results, the operator should carefully read this entire manual and fully understand all aspects of each assay step before running the assay. In addition, proper training as well as instrument maintenance is critical for obtaining optimal results in performing SMC® assays. The following notes should be reviewed and understood before the assay is set up.

- Wipe down bench and pipettes with 70% isopropanol before use.
- It is important to allow all reagents to warm to room temperature (RT), 20-25 °C.
- Use sterile filter pipette tips and reagent trays to avoid contamination.
- Pre-wet tips (aspirate and dispense within well) twice before each transfer.
- The standards prepared by serial dilution must be used within 10 minutes of preparation.

**Note:** It is recommended that the standards are prepared as the last step prior to plate setup.

- All washing must be performed with the Wash Buffer provided.
- An orbital microplate shaker for assay plate incubation (example Boekel Scientific Jitterbug™ Shaker settings #3-5) provide maximal orbital mixing without splashing liquid or causing cross-contamination.

Jitterbug™ Shaker setting #3 ~ 750 rpm

Jitterbug™ Shaker setting #4 ~ 875 rpm

Jitterbug™ Shaker setting #5 ~1000 rpm

**Note:** If using different orbital shaker, refer to recommended rpm ranges provided for each incubation step, and adjust speeds as necessary to ensure maximal orbital mixing without splashing liquid or causing cross-contamination.

- As the SMC® assay is extremely sensitive to dust particles, do not perform the assay or plate washing under direct airflow.
- Plate must also be protected from light after adding detection.
- After the assay is complete, seal the plate before reading immediately or storing temporarily at 2-8 °C. The SMCxPRO® Immunoassay System requires the use of aluminum adhesive plate seal.
- It is not recommended to store eluted products from SMC® assays overnight at 4 °C or frozen at -80 °C for later reading as performance cannot be guaranteed.
- If SMC® Read Plate has been stored at 4 °C, plate should be left at RT for 30 minutes to 1 hour on the benchtop before reading to avoid a rapid increase in temperature within SMC® Read Plate wells. Bring to RT then centrifuge the plate at 1,100 x g for 1 minute prior to reading.
- For optimal SMCxPRO® Immunoassay System performance, perform ASSIST testing daily (ideally at beginning of the day before assay is prepared).

## Precautions

Use caution when handling biological samples. Wear protective clothing and gloves.

Components of this reagent kit contain Sodium azide as a preservative. Sodium azide is a toxic and dangerous compound when combined with acids or metals. Solutions containing Sodium azide should be disposed of properly.

## Hazard Labels

| Ingredient                    | Catalogue No. | Label                                                                             |                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                      |
|-------------------------------|---------------|-----------------------------------------------------------------------------------|--------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| IFN- $\lambda$ 1 Standard     | 02-8213-00    |  | <b>Warning.</b> Harmful if swallowed, in contact with skin or if inhaled. May cause damage to organs Brain through prolonged or repeated exposure if swallowed. May cause damage to organs like respiratory tract through prolonged or repeated exposure if inhaled. Harmful to aquatic life with long lasting effects. Do not breathe dust. Wash skin thoroughly after handling. Do not eat, drink, or smoke when using this product. Use only outdoors or in a well-ventilated area. Avoid release to the environment. Wear protective gloves/ protective clothing. IF SWALLOWED: Call a POISON CENTER/ doctor if you feel unwell. Rinse mouth. IF ON SKIN: Wash with plenty of water. Call a POISON CENTER/ doctor if you feel unwell. IF INHALED: Remove person to fresh air and keep comfortable for breathing. Call a POISON CENTER/ doctor if you feel unwell. Get medical advice/ attention if you feel unwell. Wash contaminated clothing before reuse. Dispose of contents/ container to an approved waste disposal plant. |
| IFN- $\lambda$ 1 Coated Beads | 02-2213-00    | No Label Required                                                                 | Harmful to aquatic life with long lasting effects. Avoid release to the environment. Dispose of contents/ container to an approved waste disposal plant.                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                             |

| <b>Ingredient</b> | <b>Catalogue No.</b> | <b>Label</b>                                                                      |                                                                                                                                                                                                                                                                                             |
|-------------------|----------------------|-----------------------------------------------------------------------------------|---------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Assay Buffer      | 02-9950-00           | No Label Required                                                                 | Harmful to aquatic life with long lasting effects. Avoid release to the environment. Dispose of contents/ container to an approved waste disposal plant.                                                                                                                                    |
| Standard Diluent  | 02-0225-02           |  | <b>Warning.</b> May cause damage to organs Respiratory Tract through prolonged or repeated exposure if inhaled. Do not breathe dust/ fume/ gas/ mist/ vapours/ spray. Get medical advice/ attention if you feel unwell. Dispose of contents/ container to an approved waste disposal plant. |
| 10X Wash Buffer   | 02-0001-03           |  | <b>Warning.</b> Causes serious eye irritation. Harmful to aquatic life with long lasting effects. Avoid release to the environment. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing.                        |
| Elution Buffer B  | 02-0211-02           | No Label Required                                                                 | Harmful to aquatic life. Avoid release to the environment. Dispose of contents/ container to an approved waste disposal plant.                                                                                                                                                              |

# Assay Preparation

## Sample Preparation

### Preferred Method

1. Stack the filter plate on top of a 96-well receptacle plate.
2. Place 250  $\mu$ L of sample into a filter plate well and spin for  $\geq 10$  minutes at 1,100  $\times g$ .

### Optional Method

1. Centrifuge samples at  $> 13,000 \times g$  for 10 minutes immediately prior to use.
2. Carefully pipette the supernatant into a clean microcentrifuge tube, avoiding particulates and slowly aspirating below the lipid layer.

## Sample Dilution

1. Dilute the clarified samples 1:2 using the Standard Diluent (For triplicates, transfer 200  $\mu$ L of clarified sample to the sample preparation plate and add 200  $\mu$ L Standard Diluent).
2. 100  $\mu$ L per well of 1:2 diluted serum or plasma sample should be used.
3. If further sample dilution is required, samples can be diluted with the provided Standard Diluent.

## Reagent Preparation

1. Warm all reagents to RT prior to use.
2. Store the Detection Antibody away from light until ready to use.
3. Prepare 1X Wash Buffer (from 10X Wash Buffer) as follows:
  - a. Pour all 3 bottles of 10X Wash Buffer (containing 50 mL each for 150 mL total) into a container capable of holding at least 2 L. Add 1.35 L of deionized water.
  - b. Mix thoroughly by gentle inversion or with a clean, sterile stir bar.

**NOTE:** 1X Wash Buffer may be filter sterilized (refer to [Kit Storage](#)).

4. Mix IFN- $\lambda$ 1 Antibody Coated Beads on a rotisserie spin rotator, or manually by repeat inversion, for  $\geq 20$  minutes until all beads are resuspended.

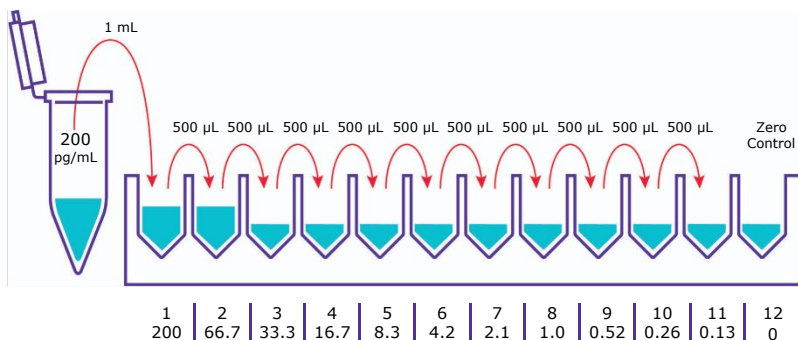
## Initial Standard Stock Preparation

1. Reconstitute lyophilized standard in 250  $\mu\text{L}$  of deionized water. Invert the vial several times to mix. Gently pulse vortex the vial for 10 seconds. Allow the vial to sit for 5-10 minutes.
2. Refer to the standard value assignment on the Certificate of Analysis for the starting concentration of the IFN- $\lambda$ 1 Standard in the vial.
3. Perform the necessary dilutions in Standard Diluent to achieve the final working concentration of 200 pg/mL in a 1.0 mL final volume.

## Standard Curve

Prepare the standard curve in a 12-channel reagent reservoir. Perform a 1:3 serial dilution of the 200 pg/mL Standard 1 for Standard 2 followed by 1:2 serial dilutions for standards 3 through 11 to achieve a curve from 200 pg/mL to 0.13 pg/mL. Standard 12 is the Blank (Standard Diluent only).

Run the standards in triplicate.



**NOTE:** Pipette gently into reservoir wells to avoid creating bubbles.

1. Add 1000  $\mu\text{L}$  Standard Diluent to well 2 of a 12-channel reservoir dilution plate
2. Add 500  $\mu\text{L}$  Standard Diluent to wells 3 through 12 of a 12-channel reservoir dilution plate.
3. Transfer 1000  $\mu\text{L}$  200 pg/mL working stock (Standard 1) into well 1.
4. Transfer 500  $\mu\text{L}$  from well 1 into well 2, mixing thoroughly. Continue serial dilutions from well 2 stopping at well 11, mixing thoroughly each time. Use a fresh tip with each transfer.

# Assay Procedure

## Target Capture

1. Pipette 100  $\mu$ L per well of Standards or 1:2 diluted Samples to assay plate.
2. Retrieve the IFN- $\lambda$ 1 Coated Bead vial from the rotator and transfer its full contents to 11.0 mL of supplied Assay Buffer. Rinse the bead vial with 0.55 mL of fresh Assay Buffer and ensure that all beads have been transferred from the original vial. Mix by gentle inversion. There should be a total volume of 12.1 mL of diluted IFN- $\lambda$ 1 Coated Beads.
3. Using a multichannel pipette, add 100  $\mu$ L per well of diluted IFN- $\lambda$ 1 Coated Beads into the assay plate.
4. Seal the assay plate with clear adhesive plate seal, applying pressure to the seal to prevent leaking and cross-contamination.
5. Incubate for 2 hours at 25 °C on microplate incubator/shaker (Jitterbug™ setting #3).
6. A minimum of 10 minutes prior to the end of target capture incubation, prepare the IFN- $\lambda$ 1 detection antibody working stock:
7. Prepare 1X detection antibody by adding 250  $\mu$ L of 20X detection antibody into 4,750  $\mu$ L of Assay Buffer and filter the diluted detection antibody using the syringe with a 0.2  $\mu$ m filter into a clean tube.
8. When incubation is complete, centrifuge the assay plate at 1,100 x g for 1 minute, place the plate on the washer magnet, and carefully remove clear adhesive plate seal to avoid splashing.

## Post-Capture Wash

Wash plate once with a plate washer using the Post Capture Wash (POSTCAP) program on the BioTek® 405 TSUVS washer.

If using automation, please contact your technical service representative for the appropriate automation procedure.

## Detection Antibody Incubation

1. After removal from the plate washer, place the assay plate onto the sphere mag plate and allow beads to form a tight pellet at the well corners for 2 minutes.
2. Using a multichannel pipette, dispense 20  $\mu$ L per well of IFN- $\lambda$ 1 Detection Antibody using reverse pipetting without disturbing the bead pellets.
3. Seal the assay plate with a new clear adhesive plate seal. Apply pressure to the seal to prevent leaking and cross-contamination.
4. Incubate for 1 hour at 25 °C on microplate incubator/shaker (Jitterbug™ setting #5). Ensure plate is protected from light during this incubation.
5. When incubation is complete, centrifuge at 1,100  $\times$  g for 1 minute then carefully remove the clear adhesive plate seal to avoid splashing.

## Post-Detection Wash

Wash the assay plate with wash buffer using the 4 cycle Pre-Transfer (4CYCPRE) program on the BioTek® 405 TSUVS washer.

If using automation, please contact your technical service representative for the appropriate automation procedure.

## Post-Detection Shake

1. After 4 cycle Pre-Transfer wash, visually verify that each well contains ~200  $\mu$ L of wash buffer.
2. Seal the assay plate with a new clear adhesive plate seal. Apply pressure to the seal to prevent leaking and cross-contamination.
3. Place the plate on the microplate/incubator shaker (Jitterbug™ setting #3) for 2 minutes. **Ensure plate is protected from light during this incubation.**
4. Remove the plate from the shaker, and centrifuge at 1,100  $\times$  g for 1 minute. Carefully remove clear adhesive plate seal to avoid splashing and place it on the plate washer to perform Final Aspiration.

## Final Aspiration

Perform the Final Aspirate program (FINASP) program on the BioTek® 405 TSUVS washer.

If using automation, please contact your technical service representative for the appropriate automation procedure.

## Elution

1. After removal from the plate washer, place the assay plate onto the sphere mag plate and allow beads to form a tight pellet at the well corners for 2 minutes.
2. Dispense 10  $\mu$ L Elution Buffer B per well using reverse pipetting without disturbing the bead pellet.
3. Seal assay plate with a new clear adhesive plate seal. Apply pressure to the seal to prevent leaking and cross-contamination.
4. Incubate the plate for 10 minutes at 25 °C on microplate incubator/shaker (Jitterbug™ setting #5). **Ensure plate is protected from light during this incubation.**
5. When incubation is complete, centrifuge at 1,100 x g for 1 minute.

## Assay Reading

### To read on the SMCxPRO® Immunoassay System

1. Place the assay plate with Elution Buffer B onto the sphere mag plate and allow beads to form a tight pellet for 2 minutes.
2. Keeping the assay plate on the magnet, carefully remove the adhesive plate seal. Using a multichannel pipette, add 10  $\mu$ L of Buffer D to center of wells containing Elution Buffer B. Use a fresh tip with each dispense.
3. Set a manual 12-channel pipette to 18  $\mu$ L and put 12 tips onto the pipettor. Transfer 18  $\mu$ L of neutralized eluate solution per well to corresponding wells of the reading plate, placed over the included plate holder by aspirating directly from the v-bottom of the plate, avoiding the pelleted beads, and changing tips with each dispensed row.
4. Seal the reading plate with new clear adhesive plate seal. Centrifuge plate for 1 minute at RT, approximately 1,100 x g. Remove the seal, inspect reading plate wells and remove bubbles if they are present.
5. Firmly seal the reading plate with aluminum plate seal using the recommend plate roller.
6. Remove the plate holder from the sealed reading plate and load it onto the SMCxPRO® Immunoassay System. Start read.

**Note:** There is a warmup period of up to 30 minutes to equilibrate plate temperature to internal instrument temperature. Once achieved the read will start automatically.

## SMC<sup>®</sup> Assay Overview

1. Prepare all reagents, standard curve, and samples as instructed.
2. Add 100  $\mu$ L of Standard/1:2 diluted samples and 100  $\mu$ L of Coated Beads to assay plate.
3. Seal and incubate for 2 hours at 25 °C on appropriate microplate incubator/shaker.



2 hours at 25 °C

4. After capture incubation, centrifuge assay plate at 1,100  $\times g$  for 1 minute.
5. Perform Post-Capture Wash.
6. Remove from washer magnet and add 20  $\mu$ L of Detection Antibody per well.
7. Seal assay plate and incubate for 1 hour at 25 °C on microplate incubator/shaker.



1 hour at 25 °C

8. Perform Post-Detection Wash.
9. Seal the assay plate and perform the post-detection shake for 2 minutes on microplate incubator/shaker.
10. Perform the Final Aspiration.
11. Remove from washer magnet and add 10  $\mu$ L of Elution Buffer B to each well of assay plate.
12. Seal assay plate and incubate for 10 minutes at 25 °C on microplate incubator/shaker.



10 minutes at 25 °C

13. Add 10  $\mu$ L of Buffer D to neutralize the eluted antibody.
14. Transfer 18  $\mu$ L of neutralized eluate to the **SMC® Read Plate**.
15. Seal **SMC® Read Plate** with aluminum adhesive plate seal for SMCxPRO® System.
16. Load on SMCxPRO® System

## Assay Characteristics

### Sensitivity

Assay sensitivity measures the true limit of quantitation of an analyte and is often defined by the Lower Limit of Quantification (LLOQ). LLOQ is calculated as the lowest concentration that can achieve CVs of < 20% and the percent recovery of the standard point is still between 80%-120%. The LLOQ of IFN- $\lambda$ 1 is 0.26 pg/mL. Please note that the published LLOQ is data generated during kit verification and can have minor variation between kit lots. For lot specific LLOQ, please see the certificate of analysis.

### Precision

The assay variations of SMC® Human IFN-  $\lambda$ 1 Immunoassay kits were studied using five normal plasma samples run in triplicate by 3 different operators on 3 different days.

- Mean intra-assay variation was 5%.
- Mean inter-assay variation was 9%.

### Cross-Reactivity

Cross-reactivity to the following analytes were tested with the following results:

- |                                        |                                         |
|----------------------------------------|-----------------------------------------|
| • IFN- $\alpha$ 2 – not cross-reactive | • IFN- $\omega$ – not cross-reactive    |
| • IFN- $\beta$ – not cross-reactive    | • IFN- $\lambda$ 2 – not cross-reactive |
| • IFN- $\epsilon$ – not cross-reactive | • IFN- $\lambda$ 3 – not cross-reactive |
| • IFN- $\gamma$ – not cross-reactive   |                                         |

Specificity to the following species samples were tested with the follow results:

- |                                                      |                               |
|------------------------------------------------------|-------------------------------|
| • Non-Human Primate (Cyno)<br>– not cross-reactive   | • Mouse – not cross-reactive  |
| • Non-Human Primate (Rhesus)<br>– not cross-reactive | • Feline – not cross-reactive |
| • Rat – not cross-reactive                           | • Canine – not cross-reactive |

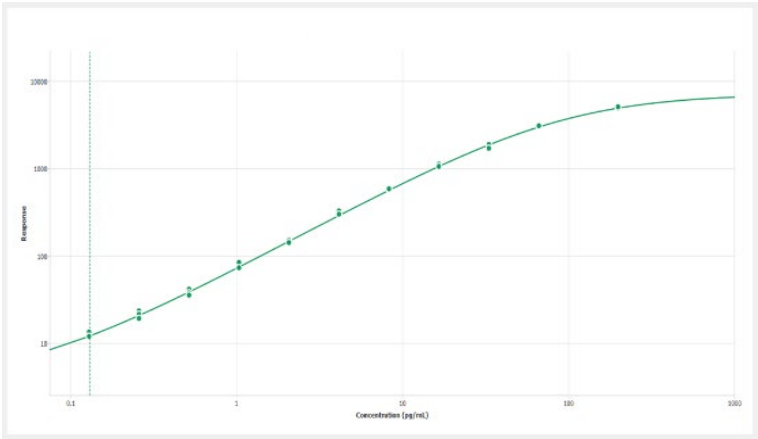
## Spike Recovery

The data represent mean percent recovery of three different concentrations of standard spiked into samples (n = 5 serum samples, 5 plasma samples).

| Sample ID | Serum Recovery | Plasma Recovery |
|-----------|----------------|-----------------|
| Sample 1  | 95%            | 103%            |
| Sample 2  | 82%            | 91%             |
| Sample 3  | 78%            | 83%             |
| Sample 4  | 70%            | 94%             |
| Sample 5  | 101%           | 102%            |
| Average   | 85%            | 95%             |

## Graph of Typical Reference Curve

Typical SMCxPRO® IFN-λ1 Immunoassay Standard Curve is shown below, not to be used to calculate data.



# Troubleshooting

| Problem                    | Probable Cause                                                           | Solution                                                                                                                                                                                                                                                                                                                                                   |
|----------------------------|--------------------------------------------------------------------------|------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Background is too high     | Background wells were contaminated                                       | <p>Avoid cross-well contamination by using seal appropriately. Pipette with multichannel pipets without touching reagent in plate. Change tips when adding reagents if cross contamination is expected.</p> <p>Ensure reagents (including wash buffer) are not contaminated.</p> <p>Insufficient washes—washer may need to be cleaned or reprogrammed.</p> |
|                            | Plate was over-incubated                                                 | Confirm plate incubation times are as recommended, particularly for the Detection incubation.                                                                                                                                                                                                                                                              |
|                            | Multichannel pipet may not be calibrated                                 | Calibrate pipets.                                                                                                                                                                                                                                                                                                                                          |
| Sample variability is high | Plate washing was not uniform                                            | Confirm that there is no residual left in the wells following post-capture wash step and Final Aspirate. Ensure that you have < 2 $\mu$ L or residual remaining in the well.                                                                                                                                                                               |
|                            | Samples may have high particulate matter or other interfering substances | Samples should be filtered according to the Assay Preparation section. Unprocessed samples could lead to higher imprecision.                                                                                                                                                                                                                               |
|                            | Plate agitation was insufficient                                         | Plate should be agitated during all incubation steps using an orbital plate shaker at a speed where beads are in constant motion without causing splashing (See <a href="#">Jitterbug™ Shaker setting</a> in Assay Best Practices section).                                                                                                                |
|                            | Cross-well contamination                                                 | Ensure that the plate is sealed well at each incubation step. If splashing occurs on plate seal, centrifuge plate at 1,100 $\times g$ for 1 minute to remove material prior to removing the seal. A new plate seal should be used every time the plate is sealed.                                                                                          |
|                            |                                                                          | Care should be taken when using same pipet tips that are used for reagent additions and that pipet tip does not touch reagent in plate.                                                                                                                                                                                                                    |

| <b>Problem</b>                                           | <b>Probable Cause</b>                                                | <b>Solution</b>                                                                                                                                                                                                                   |
|----------------------------------------------------------|----------------------------------------------------------------------|-----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Beads are lost during the wash                           | Plate washer needs optimization/cleaning                             | Contact Tech Support or local Specialist to schedule washer programming. Refer to user guide for cleaning procedure.                                                                                                              |
|                                                          | Insufficiently primed washer                                         | Washer should be primed with wash buffer prior to running the post capture wash protocol.                                                                                                                                         |
|                                                          | Beads came in contact with water                                     | Washer should be primed with wash buffer sufficiently prior to plate wash. Viscosity of water changes the performance of the magnetic particles.                                                                                  |
|                                                          | Proper magnet was not used                                           | Ensure that the SMC® magnetic plate shipped with the BioTek® 405 TSUVS Plate Washer was present on plate wash stage prior to running wash protocol.                                                                               |
| Published LLoQ was not achieved                          | Improper dilution/reconstitution of the standard reference material  | Confirm appropriate kit protocol was followed when preparing standard curve.                                                                                                                                                      |
|                                                          |                                                                      | Check plate washer to confirm no beads were lost during washes and that plate contains < 2 µL following the post-capture and final aspiration protocols.<br><br>Ensure standards are prepared before starting capture incubation. |
| Microparticles do not resuspend into homogenous solution | Beads were not properly stored and may have been frozen              | Labelled microparticles should be stored at 4 °C. If microparticles are frozen, they will not resuspend properly.                                                                                                                 |
|                                                          | Samples may be causing interference due to excess particulate matter | Samples should be properly processed prior to testing to remove particulate matter or lipids.                                                                                                                                     |

## Well Plate

|   | 1             | 2             | 3             | 4             | 5             | 6             | 7             | 8             | 9             | 10             | 11             | 12             |
|---|---------------|---------------|---------------|---------------|---------------|---------------|---------------|---------------|---------------|----------------|----------------|----------------|
| A | Standard<br>1 | Standard<br>2 | Standard<br>3 | Standard<br>4 | Standard<br>5 | Standard<br>6 | Standard<br>7 | Standard<br>8 | Standard<br>9 | Standard<br>10 | Standard<br>11 | Standard<br>12 |
| B | Standard<br>1 | Standard<br>2 | Standard<br>3 | Standard<br>4 | Standard<br>5 | Standard<br>6 | Standard<br>7 | Standard<br>8 | Standard<br>9 | Standard<br>10 | Standard<br>11 | Standard<br>12 |
| C | Standard<br>1 | Standard<br>2 | Standard<br>3 | Standard<br>4 | Standard<br>5 | Standard<br>6 | Standard<br>7 | Standard<br>8 | Standard<br>9 | Standard<br>10 | Standard<br>11 | Standard<br>12 |
| D | Sample<br>1   | Sample<br>1   | Sample<br>1   | Sample<br>2   | Sample<br>2   | Sample<br>2   | Sample<br>3   | Sample<br>3   | Sample<br>3   | Etc.           | Etc.           | Etc.           |
| E |               |               |               |               |               |               |               |               |               |                |                |                |
| F |               |               |               |               |               |               |               |               |               |                |                |                |
| G |               |               |               |               |               |               |               |               |               |                |                |                |
| H |               |               |               |               |               |               |               |               |               |                |                |                |

For research use only. Not for use in diagnostic procedures.

---

## Terms of Sale

THIS PRODUCT IS INTENDED FOR USE BY AN ACADEMIC OR NOT-FOR-PROFIT INSTITUTION TO BE USED FOR ACADEMIC AND/OR NOT-FOR-PROFIT RESEARCH, WHICH IS FURTHER DEFINED BELOW. FOR COMMERCIAL USE PLEASE CONTACT US AT THE E-MAIL ADDRESS BELOW. BY OPENING THIS PRODUCT, YOU ("PURCHASER") HEREBY REPRESENT THAT YOU HAVE THE RIGHT AND AUTHORITY TO LEGALLY BIND YOURSELF AND/OR YOUR EMPLOYER INSTITUTION, AS APPLICABLE, AND CONSENT TO BE LEGALLY BOUND BY THE TERMS OF THIS ACADEMIC USE AGREEMENT. IF YOU DO NOT AGREE TO COMPLY WITH THESE TERMS, YOU MAY NOT OPEN OR USE THE PRODUCT AND YOU MUST CALL CUSTOMER SERVICE (1-800-645-5476) TO ARRANGE TO RETURN THE PRODUCT FOR A REFUND.

"PRODUCT" means SMCxPRO® Immunoassay Instrument (95-0100-00, 70-0100-00, 95-0100-00-JPN).

"Commercial Product" means any product intended for: (i) current or future sale; (ii) use in a fee-for-service; or (iii) any diagnostic, clinical, or therapeutic use.

"Academic or Not-For-Profit Research" means any internal in vitro research use by individuals employed by an academic or not-for-profit institution. Such research specifically excludes the following uses of whatever kind or nature:

- Re-engineering or copying the PRODUCT
- Making derivatives, modifications, or functional equivalents of the PRODUCT
- Obtaining patents or other intellectual property rights claiming use of the PRODUCT
- Using the PRODUCT in the development, testing, or manufacture of a Commercial Product
- Using the PRODUCT as a component of a Commercial Product
- Reselling or licensing the PRODUCT
- Using the PRODUCT in clinical or therapeutic applications including producing materials for clinical trials
- Using the PRODUCT to provide a service to any third party
- Using the PRODUCT in collaboration or to enable a commercial entity
- Commercial Use of the PRODUCT to make a "home brew" assays, which includes a LDT (Lab Developed Test) assay(s) or any related commercial testing
- Transfer of PRODUCT to a site that is not the site where the Product was originally installed.

---

Access to the PRODUCT is limited solely to those officers, employees and students of PURCHASER's not-for-profit institution who need access to the PRODUCT for internal in vitro research use. PURCHASER shall comply with all applicable laws in its use and handling of the PRODUCT and shall keep it under reasonably safe and secure conditions to prevent unauthorized use or access.

These use restrictions will remain in effect for as long as PURCHASER possesses the PRODUCT.

COMMERCIAL OR NON-ACADEMIC ENTITIES INTERESTED IN PURCHASING OR USING THE PRODUCT MUST CONTACT [limitedcommercial@milliporesigma.com](mailto:limitedcommercial@milliporesigma.com) AND AGREE TO SEPARATE TERMS OF LIMITED COMMERCIAL USE PRIOR TO USE OR PURCHASE.

## Notice

We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

The information in this document is subject to change without notice and should not be construed as a commitment by the manufacturing or selling entity, or an affiliate. We assume no responsibility for any errors that may appear in this document.

## Technical Assistance

Visit the tech service page at [SigmaAldrich.com/techservice](https://SigmaAldrich.com/techservice).

## Terms and Conditions of Sale

Warranty, use restrictions, and other conditions of sale may be found at [SigmaAldrich.com/terms](https://SigmaAldrich.com/terms).

## Safety Data Sheets (SDS)

Safety Data Sheets are available on the product page at [SigmaAldrich.com](https://SigmaAldrich.com).

## Contact Information

For the location of the office nearest you, go to [SigmaAldrich.com/offices](https://SigmaAldrich.com/offices).

The life science business of Merck operates as MilliporeSigma in the U.S. and Canada.

Merck, Millipore, Stericup, SMCxPRO, Millex, Multiscreen and Sigma-Aldrich are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.

© 2024 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved.

For research use only. Not for use in diagnostic procedures.