

SMC™ Human IL-17F High Sensitivity Immunoassay Kit

Microparticle Assay

Catalog # 03-0164-00

SMC™ Human IL-17F High Sensitivity
Immunoassay Kit for the Quantitative Determination
of IL-17F in Human Plasma and Serum

FOR RESEARCH USE ONLY

NOT FOR USE IN DIAGNOSTIC PROCEDURES

Manufactured & Distributed by:



EMD Millipore Corporation
3050 Spruce Street
St. Louis, Missouri 63103
United States of America
emdmillipore.com

The M logo is a trademark of Merck KGaA, Darmstadt, Germany.
© 2013 EMD Millipore Corporation,
Billerica, MA 01821 USA.

This page intentionally left blank.

TABLE OF CONTENTS

INTRODUCTION	2
SUPPLIES	3
Reagents Provided	3
Storage Instructions	3
Required Supplies Not Provided	4
TECHNICAL HINTS	7
PRECAUTIONS	8
ASSAY PREPARATION	9
Reagent Preparation	9
Sample Preparation	9
Initial Standard Stock Preparation	10
Standard Curve	10
ASSAY PROCEDURE	11
Target Capture	11
Post-Capture Wash	11
Detection	12
Post Detection Wash	12
Post Detection Shake	12
Final Aspiration	12
Elution	12
ASSAY READING	13
To Read On Erenna® Immunoassay System	13
To Read On SMCxPRO™ Immunoassay System	13
APPENDIX A: SMC™ QUICK ASSAY GUIDE	14
GRAPH OF TYPICAL REFERENCE CURVE	15
ASSAY CHARACTERISTICS	16
TROUBLESHOOTING GUIDE	17
ORDERING INFORMATION	19
WELL MAP	20

INTRODUCTION

The Single Molecule Counting (SMC™) Human IL-17F High Sensitivity Immunoassay uses a quantitative fluorescent sandwich immunoassay technique to measure IL-17F in human plasma and serum samples. A capture antibody specific for human IL-17F has been pre-coated onto paramagnetic microparticles (beads). The user pipettes beads, standards, and samples into uncoated microplate wells. During incubation, the IL-17F 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 IL-17F that has been captured onto the beads, thus completing the immunosandwich. Elution buffer is then added and incubated. The elution buffer dissociates the bound protein sandwich from the beads surface releasing the labeled antibodies. These antibodies are separated during transfer to a final microplate. The plate is loaded into the Erenna® or SMCxPRO™ System where the labeled molecules are detected and counted. The number of fluor-labeled detection antibodies counted is directly proportional to the amount of IL-17F present in the sample when captured. The amount of IL-17F in unknown samples is interpolated from a standard curve.

**FOR RESEARCH USE ONLY
NOT FOR USE IN DIAGNOSTIC PROCEDURES**

SUPPLIES

The SMC™ Human IL-17F High Sensitivity Immunoassay Kit includes all reagents listed in *Table 1: Reagents Provided*. Additional reagents and supplies are required to run this immunoassay and are listed in *Table 2: Additional Supplies Required (not provided)*. All reagents supplied are for Research Use Only.

Table 1: Reagents Provided

Item #	Description	Shipping Conditions	Storage Conditions	Component Part No.	Packaging Details
1	Assay Buffer	With cold pack	2 - 8°C	02-9994-00	2 x 20 mL
2	Human IL-17F Coated Beads	With cold pack	2 - 8°C	02-2164-00	1 x 550 µL
3	Standard Diluent	With cold pack	2 - 8°C	02-0225-02	2 x 20 mL
4	Human IL-17F Detection Antibody	With cold pack	2 - 8°C	02-1164-00	1 x 270 µL
5	Human IL-17F Standard	With cold pack	2 - 8°C	02-8164-00	1 lyophilized vial
6	10X Wash Buffer	With cold pack	2 - 8°C	02-0001-03	1 x 50 mL
7	Buffer D	With cold pack	2 - 8°C	02-0446-00	1 x 6 mL
8	Elution Buffer B	With cold pack	2 - 8°C	02-0211-02	1 x 5 mL

Storage Instructions

The SMC™ Human IL-17F 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, EMD Millipore PN SCGPU11RE 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.

Table 2: Additional Supplies Required (not provided)**Instrumentation**

Item #	Product Description	Supplier	Product Number	Product Uses
1	12-Channel Manual Pipette 10 – 20 µL	--	--	Transferring 10 µL
2	12-Channel Manual Pipette 20 – 250 µL	--	--	Transferring 20 µL, 100 µL
3	Tube Rotator	--	--	Microparticle resuspension
4	Sphere Mag Plate	EMD Millipore	90-0003-02	Capturing/pelleting microparticles
5	Jitterbug™ Microplate Incubator/Shaker	EMD Millipore	70-0009-00	Incubating/Shaking at 25°C
6	VWR® Microplate Shaker	VWR International	12620-926	Plate shaking for overnight incubation, if recommended
7	Bio-Tek ELx™ 405 Microplate Washer	EMD Millipore	95-0004-05	Automated plate washing option
8	Tecan Hydroflex™ Microplate Washer	EMD Millipore	95-0005-02	Automated plate washing option
9	Centrifuge able to reach speed of 1,100 x g	--	--	Centrifuging samples, plates
10	Micro-Centrifuge	--	--	Centrifuge samples, provided Detection Antibody
11	ALPS™ 50V Microplate Heat Sealer	EMD Millipore	70-0018-00	Heat sealing 384- well plates before Erenna Reading

Additional Supplies Required (not provided) continued

Materials

Item #	Product Description	Supplier	Product Number	Product Uses
12	12-Channel Reagent Reservoir (sterile)	Argos/Cole Parmer	04395-33	Standard curve dilution
13	VistaLab™ 25 mL Reservoirs	Fisher Scientific	21-381-27C	Addition of Reagents
14	MultiScreenHTS BV 96-Well Filter Plate	EMD Millipore	MSBVN1210	Sample filtration
15	96-well V-bottom plate	Fisher Scientific	14-222-241	Assay plate
16	5 mL Luer-Lok™ Syringe	Fisher Scientific	14-829-45	Detection Antibody filtration
17	0.2 µm Syringe Filter	EMD Millipore	SLGPR33RS	Detection Antibody filtration
18	Nunc™ Clear Adhesive Plate Seal	Fisher Scientific	236366	Sealing assay plate
19	384-well round bottom plates	Fisher Scientific	12-565-384	Erenna® reading plate
20	Heat sealing foil	Fisher Scientific	NC0276513	Sealing plates for Erenna® reading
21	1L Stericup® Filter; 0.22 µm	EMD Millipore	S2GPU11RE	Filter sterilizing Erenna® system buffer
22	SMCxPRO™ 384-well plate, 1 plate with adhesive seal	EMD Millipore	02-1008-00	SMCxPRO™ reading plate, seal
23	SMCxPRO™ 384-well plate, case of 32	EMD Millipore	ABB2-00160A	SMCxPRO™ reading plate
24	SMCxPRO™ aluminum adhesive plate seals	Fisher Scientific	276014	SMCxPRO™ reading plate seals
25	Plate Roller	Fisher Scientific	NC9185793	Creates secure/even seal for each well of SMCxPRO™ reading plate

Additional Supplies Required (not provided) continued

Reagents

Item #	Product Description	Supplier	Product Number	Product Uses
26	Universal plate cover	Fisher Scientific	253623	Covers assay plate
27	500 mL Container	--	--	Wash Buffer Dilution
28	Micro-centrifuge tubes	--	--	Sample storage, standard preparation
29	Elution Buffer (5 mL)	EMD Millipore	02-0002-04	Required for Erenna® maintenance
30	SMC™ 10X Wash Buffer (1 L)	EMD Millipore	02-0111-00	Automated plate washing
31	SMC™ 10X System/wash Buffer with Proclin (1 L)	EMD Millipore	02-0111-03	Use in Erenna® platform
32	De-ionized or Distilled water	--	--	Dilution of 10X Wash or System Buffer

Please contact your technical services representative for additional information or assistance selecting required but not provided supplies.

TECHNICAL HINTS

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. The following notes should be reviewed and understood before the assay is set up.

Assay Hints

1. Wipe down bench and pipettes with 70% isopropanol before use.
2. It is important to allow all reagents to warm to room temperature (20 - 25°C).
3. Use sterile filter pipette tips and reagent trays to avoid contamination.
4. Pre-wet tips (aspirate and dispense within well) twice before each transfer.
5. The standards prepared by serial dilution must be used within 10 minutes of preparation.
 - a. It is recommended that the standards are prepared as the last step prior to plate setup.
6. All washing must be performed with the wash buffer provided.
7. The recommended plate shaker settings are between #3 - #7 to provide maximal orbital mixing without splashing liquid or causing cross-contamination.
8. After the assay is complete, the plate should be read immediately.
 - a. For Erenna® Immunoassay System, use heat sealing plate foil.
 - b. For SMCxPRO™ Immunoassay System – use adhesive seal.
9. The plates may be stored at 2-8°C for up to 48 hours away from light if same day reading is not possible.
 - a. After the assay is complete, seal the plate before storing at 2 - 8°C
 - i. For Erenna® Immunoassay System, use heat sealing plate foil
 - ii. For SMCxPRO™ Immunoassay System – use aluminum adhesive plate seal
 - b. Bring to RT then centrifuge the plate at 1,100 $x g$ for 1 minute prior to reading.

Instrument Hints





10. For optimal Erenna® performance, execute the following prime of the instrument before reading:
 - a. Cycle routine (10,000 μL at 1,000 $\mu L/min$)
 - b. Bubble test (200 μL at 1,000 $\mu L/min$)
 - c. Complete Erenna® calibration prior to reading the plate.

Note: If carry-over is experienced: perform a clean routine using a 384-well plate and 20 $\mu L/well$:

- i. 3 wells of elution buffer
 - ii. 1 well of 10% bleach
 - iii. 5 wells of elution buffer
11. For optimal SMCxPRO™ performance, perform ASSIST testing on a daily basis (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 approximately 0.08% 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.

Ingredient, Cat #		Full Label	
10X Wash Buffer	02-0001-03		<p>Warning. 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.</p>
Human IL-17F Standard	02-8164-00	 	<p>Warning. Harmful if swallowed. May cause damage to organs Respiratory Tract through prolonged or repeated exposure if inhaled. Do not breathe dust/ fume/ gas/ mist/ vapors/ spray. Wash skin thoroughly after handling. Do not eat, drink or smoke when using this product. IF SWALLOWED: Call a POISON CENTER/doctor if you feel unwell. Get medical advice/ attention if you feel unwell. Rinse mouth. Dispose of contents/ container to an approved waste disposal plant.</p>
Human IL-17F Beads	02-2164-00	No symbol required.	Harmful to aquatic life with long lasting effects. Avoid release to the environment.
Standard Diluent	02-0225-02		<p>Warning. May cause damage to organs Respiratory Tract through prolonged or repeated exposure if inhaled. Do not breathe dust/ fume/ gas/ mist/ vapors/ spray. Get medical advice/ attention if you feel unwell. Dispose of contents/ container to an approved waste disposal plant.</p>

ASSAY PREPARATION

Reagent Preparation

1. Warm all reagents to room temperature (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 50 mL of 10X Wash Buffer into a container capable of holding at least 500 mL. Add 450 mL 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 Storage Instructions*)

4. Mix IL-17F Antibody Coated Beads on a rotisserie spin rotator, or manually by repeat inversion, for ≥ 20 minutes until all beads are resuspended.

Sample Preparation

1. Prepare samples by one of the following methods:
 - If using a microcentrifuge: Centrifuge samples at $>13,000 \times g$ for 10 minutes immediately prior to use. Carefully pipette the supernatant into a clean microcentrifuge tube, avoiding particulates and slowly aspirating below the lipid layer.
 - If using a filter plate with prefilter: Stack the filter plate on top of a 96-well receptacle plate. Place 250 μ L of sample into a filter plate well and spin for ≥ 10 minutes at $1,100 \times g$.
2. Sample dilution:
 - Dilute the clarified samples 1:2 using the Standard Diluent (*e.g. for triplicates, transfer 250 μ L of clarified sample to the sample preparation plate and add 250 μ L Standard Diluent*).
 - 100 μ L per well of 1:2 diluted plasma or serum should be used.

ASSAY PREPARATION (continued)

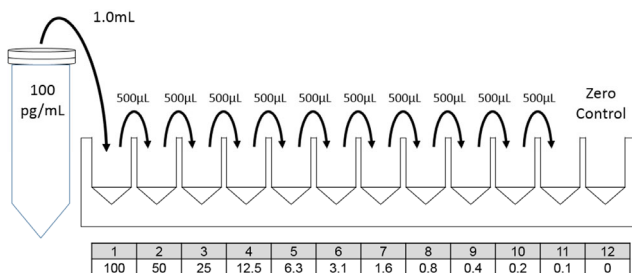
Initial Standard Stock Preparation

1. Reconstitute lyophilized standard in 250 μL 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 IL-17F Standard in the vial.
3. Perform the necessary dilutions in Standard Diluent to achieve the final working concentration of 100 pg/mL in a 1.0 mL final volume.

Standard Curve

Prepare the standard curve in a 12-channel reagent reservoir. Perform 1:2 serial dilutions of the 100 pg/mL Standard 1 for Standards 2 through 11 to achieve a curve from 100 pg/mL to 0.1 pg/mL. Standard 12 is the Blank (*Standard Diluent only*).

Run the standards in triplicate.



1. Add 500 μL Standard Diluent to wells 2 through 12 of a 12-channel reservoir dilution plate.
2. Transfer 1,000 μL 100 pg/mL working stock (Standard 1) into well 1.
3. Transfer 500 μL 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 μ L per well of Standards or 1:2 diluted Samples to assay plate.
2. Following mixing of the coated beads, immediately before adding to the assay plate, add the entire vial of coated Beads to 11.0 mL of supplied Assay Buffer. Rinse bead vial with 0.55 mL of Assay Buffer and ensure that all beads have been transferred. Mix by gentle inversion. There should be a total volume of 12.1 mL of diluted Coated Beads.
3. Pipette 100 μ L per well of the Coated Beads into assay plate.
4. Seal assay plate with clear adhesive plate seal, apply pressure to seal to prevent leaking and cross-contamination.
5. Incubate for 2 hours at 25°C on microplate incubator/shaker (Jitterbug setting #4).

Optional

For convenience, overnight incubation (16-18 hours) at 4°C can be performed on VWR® microplate shaker (setting #5). The next day, following overnight incubation, the plate should be equilibrated to room temperature by shaking for 30 minutes on microplate incubator/shaker (Jitterbug setting #4)

6. Approximately 10 minutes prior to the end of target capture incubation, prepare the detection antibody using one of the following methods:
 - a. Centrifuge 20X detection antibody at 14,000 x g for 5 minutes. Prepare 1X detection antibody by adding 250 μ L of the centrifuged supernatant into 4,750 μ L of Assay Buffer.
 - b. Prepare 1X detection antibody by adding 250 μ L of 20X detection antibody into 4,750 μ L of Assay Buffer and filter the diluted detection antibody using the syringe with a 0.2 μ m filter into a clean tube.
7. When incubation is complete, centrifuge at 1,100 x g for 1 minute and carefully remove clear adhesive plate seal to avoid splashing.

Post-Capture Wash

Wash plate once with a plate washer.

Plate Washer

- a. BioTek; Post Capture Wash (POSTCAP) or
- b. HydroFlex; Post Capture Wash (PCW)

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

ASSAY PROCEDURE (continued)

Detection

1. After removal from plate washer, dispense 20 μ L per well of Detection Antibody without disturbing the bead pellet. *(It is recommended to change tips)*
2. Seal assay plate with clear adhesive plate seal.
3. Incubate for 1 hour at 25°C on microplate incubator/shaker (Jitterbug setting #5).
4. After incubation, carefully remove clear adhesive plate seal to avoid splashing.

Post-Detection Wash

Wash assay plate 4 times with a plate washer.

Plate Washer

- a. BioTek; 4 cycle Pre-Transfer (4CYCPRE) or
- b. HydroFlex; 4 cycle Pre-Transfer (4cyPrTra)

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 μ L of wash buffer.
2. Seal assay plate with clear adhesive plate seal and apply pressure to the seal to prevent leaking and cross-contamination.
3. Place plate on microplate/incubator shaker for 1.5 minutes (Jitterbug setting #3)
4. Remove the plate from the Jitterbug, carefully remove clear adhesive plate seal to avoid splashing and place it on the plate washer to perform Final Aspiration.

Final Aspiration

Plate Washer

- a. BioTek; Final Aspirate (FINASP)
- b. HydroFlex; Final Aspirate (FA_V1)

Elution

1. Dispense 10 μ L Elution Buffer B per well using reverse pipetting without disturbing the bead pellet. *(It is recommended to change tips)*
2. Seal assay plate with a clear adhesive plate seal
3. Incubate plate for 10 minutes at 25°C on microplate incubator/shaker (Jitterbug setting #5).

ASSAY READING

To read on the Erenna® Immunoassay System

1. Add 10 μ L per well of Buffer D using reverse pipetting to Erenna® reading plate (Fisher Scientific PN 12-565-384) using a 12-channel manual P20.
2. Place assay plate onto sphere mag plate and allow beads to form a tight pellet for ≥ 2 minutes.
3. While keeping assay plate on the sphere mag plate, gently remove clear adhesive plate seal and transfer 10 μ L of eluate from assay plate to reading plate by aspirating directly from the v-bottom of the plate, avoiding the pelleted beads, and changing tips with each dispensed row.
4. Seal reading plate with clear adhesive plate seal. Centrifuge plate for 1 minute at RT, approximately 1,100 x g.
5. Seal reading plate with heat sealing foil (Fisher Scientific PN NC0276513) according to manufacturer's instructions for the heat sealer.
6. Load completed reading plate onto the Erenna® Immunoassay System.

To read on the SMCxPRO™ Immunoassay System:

1. Add 10 μ L per well of Buffer D using reverse pipetting to a fresh 96 well assay plate, using a 12-channel manual P20.
2. Place assay plate with Elution Buffer B onto sphere mag plate and allow beads to form a tight pellet for 2 minutes.
3. While keeping the assay plate containing eluate on sphere mag plate, gently remove clear adhesive seal and transfer 10 μ L of eluate to the assay plate containing Buffer D by aspirating directly from the V-bottom of the plate, avoiding the pelleted beads, and changing tips with each dispensed row.
4. Seal this plate with a clear adhesive plate seal.
5. Place the plate (containing eluted, neutralized antibody solution) into Jitterbug and shake for 2 minutes at 25°C (Jitterbug setting #5), centrifuge plate for 1 minute at RT, approximately 1,100 X g.
6. Gently remove clear adhesive plate seal and transfer 20 μ L of neutralized eluate solution per well to corresponding wells of the SMCxPRO™ reading plate (EMD Millipore PN 02-1008-00), placed over the included plate holder.
7. Seal reading plate with clear adhesive plate seal. Centrifuge plate for 1 minute at RT, approximately 1,100 x g. Remove plate sealer, inspect reading plate wells and remove bubbles if they are present.
8. Firmly seal reading plate with aluminum adhesive plate seal using the recommended plate roller.
9. Remove the plate holder from the sealed reading plate and load it onto the SMCxPRO™ Immunoassay System. Start read.

Note: there is a smart warm up period of up to 30 minutes to wait for the read plate to be close to the internal instrument temperature. Once achieved the read will start automatically.

APPENDIX A: SMC™ Quick Assay Guide

1. Prepare all reagents, standard curve, and samples as instructed.
2. Add 100 μ L of Standard/1:2 samples and 100 μ L of **Coated Beads** to **assay plate**.
3. Seal and incubate for 2 hours at 25°C or overnight (16-18 hours) at 4°C on appropriate microplate incubator/shaker.



2 hours 25°C/overnight 4°C

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



1 hour 25°C

8. Perform Post-Detection Wash.
9. Perform Post-Detection Shake for 1.5 minutes on Jitterbug setting #3.
10. Perform Final Aspiration.
11. Remove from washer magnet and add 10 μ L of **Elution Buffer B** to each well
12. Seal and incubate for 10 minutes at 25°C on microplate incubator/shaker.



10 minutes 25°C

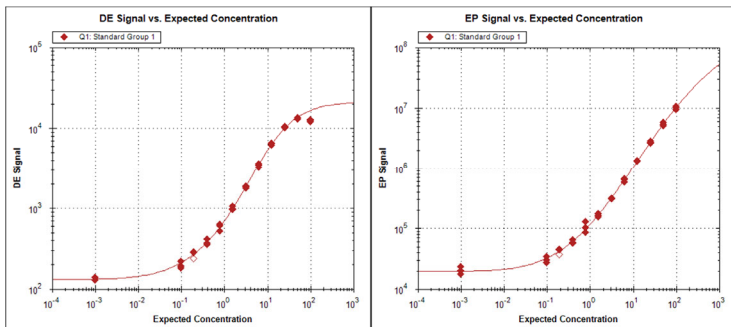
13. Neutralize eluted antibody.
14. Seal **reading plate** with pierceable foil for Erenna® or aluminum adhesive plate seal for SMCxPRO™.



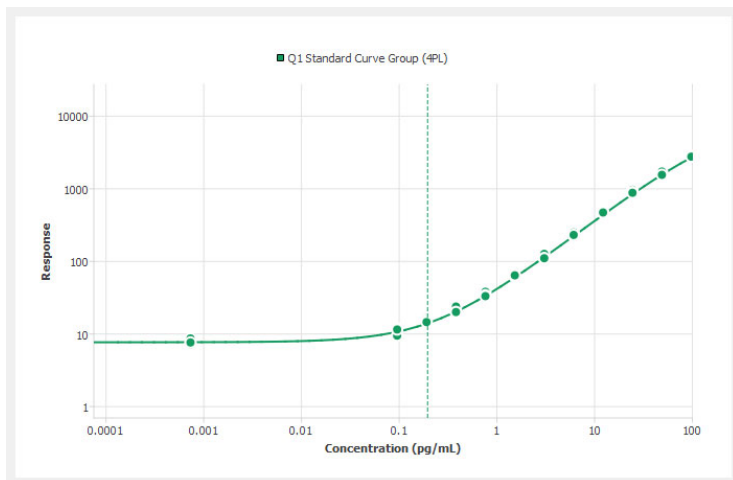
LOAD ON ERENNA® or SMCxPRO™ SYSTEM

GRAPH OF TYPICAL REFERENCE CURVE

Typical Erenna[®] Immunoassay System Standard Curve in DE and EP signal, not to be used to calculate data.



Typical SMCxPRO[™] Immunoassay System Standard Curve, not to be used to calculate data.



ASSAY CHARACTERISTICS

A. 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 IL-17F is 0.2 pg/mL. The reported value is the average of multiple assays (n = 10).

B. Precision

The assay variations of EMD Millipore's SMC™ Human IL-17F High Sensitivity Immunoassay kit were studied using three different concentrations of IL-17F (generated from serial dilutions of the SMC™ IL-17F standard).

The mean intra-assay variation (n = 20) was calculated from a single assay for each of the three concentrations of IL-17F standard. Mean intra-assay variation was < 10%.

The mean inter-assay variation (n = 5) was calculated from six independent assays of the same three dilutions of IL-17F used to determine intra-assay variation. Mean inter-assay variation was < 15%.

C. Specificity/Cross-Reactivity

Non-Human Primate samples found to be detectable

D. Spike Recovery

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

Sample ID	Recovery %
Plasma Sample 1	92
Plasma Sample 2	108
Plasma Sample 3	106
Plasma Sample 4	100
Plasma Sample 5	117
Plasma Average	105

Sample ID	Recovery %
Serum Sample 1	92
Serum Sample 2	95
Serum Sample 3	100
Serum Sample 4	97
Serum Sample 5	86
Serum Average	94

TROUBLESHOOTING GUIDE

Problem	Probable Cause	Solution
Background is too high	Background wells were contaminated	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.
		Ensure reagents (including wash and system buffers) are not contaminated.
		Insufficient washes—washer may need to be cleaned or reprogrammed.
	Instrument needs cleaning	See Technical Hints for appropriate Erenna® cleaning protocol.
	Plate was over-incubated	Confirm plate incubation times are as recommended, particularly for the Detection incubation.
Sample variability is high	Multichannel pipet may not be calibrated	Calibrate pipets.
	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 µL or residual remaining in the well.
	Samples may have high particulate matter or other interfering substances	Samples should be centrifuged or 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 a vertical plate shaker at a speed where beads are in constant motion without causing splashing (~650 - 1000 RPM).
	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 x 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.
Beads are lost during the wash	Plate washer needs optimization/cleaning	Contact Tech Support or local BCS 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.

TROUBLESHOOTING GUIDE (continued)

Problem	Probable Cause	Solution
Beads are lost during the wash (continued)	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 mag plate (EMD Millipore PN 90-0003-02) 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 μ L following the post-capture and final aspiration protocols.
		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.

ORDERING INFORMATION

To place an order or to obtain additional information about SMC™ products, please contact your Customer Service or Technical Support Specialist.

Contact information for each region can be found on our website:

emdmillipore.com/contact

Conditions of Sale

For Research Use Only. Not for Use in Diagnostic Procedures.

Safety Data Sheets (SDS)

Safety Data Sheets for EMD Millipore products may be ordered by fax or phone or through our website at emdmillipore.com/msds

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

This page intentionally left blank.



EMD Millipore Corporation
3050 Spruce Street
St. Louis, Missouri 63103
United States of America

Toll-Free US: (800) 645-5476

Fax: (800) 645-5439

emdmillipore.com