

Technical Bulletin

Duolink® flowPLA Detection Kit - Far Red

DUO94004

Storage Temperature -20 °C

Product Description

Duolink® flowPLA Detection Kit – Far Red contains all the necessary Duolink PLA® reagents to perform the amplification and detection of bound PLA probes by flow cytometry. The detection oligonucleotides contain a fluorophore ($\lambda_{\text{ex}} = 644 \text{ nm}/\lambda_{\text{em}} = 669 \text{ nm}$).

Experiments conducted using Duolink® flowPLA reagents can detect protein interactions, protein expression levels, and post-translational modifications at the single molecule level in fixed, suspended cells.

Components

Sufficient components are provided for 40 tests, based on 100 μL total reaction volume covering 100,000 cells.

- 5x Ligation Buffer
Contains oligonucleotides that hybridize to the PLA probes and all components needed for ligation except the ligase. Cat. No. DUO82009-40 TST, 800 μL
- Ligase (1 unit/ μL) Cat. No. DUO82027, 100 μL
- Polymerase (10 units/ μL) Cat. No. DUO82028, 50 μL
- 5x Amplification Buffer – Contains all components needed for rolling-circle amplification (RCA) except the polymerase. Cat. No. DUO82050-40 TST, 800 μL
- 5x flowPLA Detection Solution Far Red
Contains oligonucleotides labeled with a fluorophore that hybridize to the RCA product. Cat. No. DUO84024-40 TST, 800 μL

Reagents and Equipment Required

(But Not Provided)

To perform a complete Duolink® flowPLA experiment, one will need two primary antibodies (IHC or ICC/IF validated) that recognize two target epitopes. Additional reagents include a pair of PLA probes (one 100RXN PLUS and one 100RXN MINUS) and flowPLA detection reagents of choice. Recommended reagents include Duolink® Wash Buffers and PBS.

Precautions and Disclaimer

For R&D use only. Not for drug, household, or other uses. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

Preparation Instructions

Thaw the 5x Ligation, 5x Amplification, and 5x flowPLA Detection Far Red buffers at room temperature and vortex before use. Dilute the required volumes of each 5x solution 5-fold with ultrapure water immediately before use. Do not store diluted reagents.

Note: The 5x Ligation Buffer contains DTT that may precipitate at -20 °C. Make sure the DTT is completely dissolved and vortexed before use.

The Duolink® Detection solutions are light-sensitive. Protect from light.

The ligase and polymerase enzymes should be kept cold (-20 °C) at all times; use a freezing block when removing them from the freezer. Quick spin the vial before pipetting. Add the enzyme to the appropriate reaction mix immediately before use. Vortex the mix after addition of enzyme. Do not store diluted reagents.

Storage/Stability

Store the flowPLA reaction components at -20 °C. The enzymes should be kept cold (-20 °C) at all times, use a freezing block when removing them from the freezer.

Procedure

Note: Duolink® PLA reagent volumes are based on a 40 µL reaction volume for a 1 cm² sample on a microscope slide or a 100 µL reaction volume at ~ 1,000 cells/µL for flow cytometry. However, volumes may need to be adjusted according to the sample size or number of cells of the sample.

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.

Standard Warranty

The applicable warranty for the products listed in this publication may be found at SigmaAldrich.com/terms.

Contact Information

For the location of the office nearest you, go to SigmaAldrich.com/offices.

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

Merck 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.

© 2022 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved.
DUO94004bul Rev 05/22

