

Product Information



Prostaglandin D₂ FPIA Reagent - Red

Item No. 10010445

Laboratory Procedures

This vial contains prostaglandin D₂ (PGD₂) monoclonal antibody and a rhodamine conjugated tracer in the presence of a proprietary formulation of stabilizers. For long term storage, we suggest that the reagent be stored as supplied at -20°C; it will be stable for at least one year. Reconstitute a 400 dtn vial of PGD₂ FPIA Reagent with 12 ml of FPIA Buffer or a 2,000 dtn vial with 60 ml of FPIA Buffer. *NOTE: 12 ml of reconstituted reagent is enough for either a standard 96- or 384-well plate.* Allow the reagent to equilibrate to room temperature prior to use. Store unused PGD₂ FPIA Reagent at 4°C and use within 72 hours. For your convenience, we have supplied a 20% surplus of reagent.

FPIA Buffer

Buffer for use in the FPIA can be purchased from Cayman as a 10X concentrated stock solution (Item No. 400501).

Suggested Assay Protocol

This product has been tested and formulated to work exclusively with the PGD₂ FPIA Assay Kit from Cayman (Item No. 10007385). This product may not perform as described if used with other assay reagents or protocols.

The PGD₂ standard for this assay should be prepared in FPIA Buffer using 3-fold serial dilutions beginning at 300 ng/ml. The volumes given below are for a 96-well format; decrease the volume to 25 µl for a 384-well format. *NOTE: This is an abbreviated protocol. If you are not familiar with this assay, please contact us for a complete protocol.*

1. Add 100 µl of standard or sample to appropriate wells.
2. Add 100 µl of FPIA Reagent to each well.
3. Incubate for 60-90 minutes at room temperature.
4. Read the plate(s) with excitation and emission wavelengths of 560 nm and 595 nm, respectively. The plate reader used at Cayman also employs a dichroic filter at 505 nm. Some instruments may not utilize this type of filter. The excitation polarizer is set in the horizontal position, and the emission polarizer is dynamic. The measurements are taken with the z-height set to the middle of the well and the G-factor set to one.

Related Product

Prostaglandin D₂ FPIA Kit - Red - Item No. 10007835

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog.**

Copyright Cayman Chemical Company, 10/26/2010

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com