

PRODUCT INFORMATION



Primary Eicosanoid HPLC Mixture

Item No. 10001

Purity: ≥98% for each compound
Supplied as: A solution in methyl acetate (100 µg of each compound)
Storage: -20°C
Stability: ≥6 months

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

This mixture contains the primary COX products produced by most mammalian tissues. Contents: Prostaglandin D₂, Prostaglandin E₂, 6-keto Prostaglandin F_{1α}, Prostaglandin F_{2α}, and Thromboxane B₂ (100 µg each).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/15/2020

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM