

PRODUCT INFORMATION



Prostaglandin HPLC Mixture

Item No. 10002

Synonyms: PG HPLC Mixture
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.

Laboratory Procedures

This mixture contains primary prostaglandins produced from arachidonic acid and dihomo-γ-linolenic acid. Contents: Prostaglandin E1, Prostaglandin E2, Prostaglandin F1α, 6-keto Prostaglandin F1α, and Prostaglandin F2α (100 µg each).

Name:	CAS No.:
Prostaglandin E1	745-65-3
Prostaglandin E2	363-24-6
Prostaglandin F1α	745-62-0
6-keto Prostaglandin F1α	58962-34-8
Prostaglandin F2α	551-11-1

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, 08/05/2022

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