

PRODUCT INFORMATION



Cyclopentenone Prostaglandin HPLC Mixture

Item No. 10000

Synonym: Cyclopentenone PG HPLC Mixture
Purity: ≥98% for each compound
Supplied as: A solution in methyl acetate (100 µg of each compound)
Storage: -80°C
Stability: ≥6 months

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

Laboratory Procedures

The cyclopentenone prostaglandin HPLC mixture contains all of the major UV-absorbing cyclopentenone prostaglandins and their precursors supplied in methyl acetate. This HPLC mixture contains 100 µg of the following cyclopentane prostaglandins and precursors: prostaglandin A₂, prostaglandin B₂, prostaglandin D₂, prostaglandin E₂, prostaglandin J₂, and 15-deoxy-Δ^{12,14}-prostaglandin J₂.

Name:	CAS No.:
Prostaglandin A ₂	13345-50-1
Prostaglandin B ₂	13367-85-6
Prostaglandin D ₂	41598-07-6
Prostaglandin E ₂	363-24-6
Prostaglandin J ₂	60203-57-8
15-deoxy-Δ ^{12,14} -Prostaglandin J ₂	87893-55-8

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, 07/25/2022

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