

PRODUCT INFORMATION



Prostaglandin Metabolite HPLC Mixture

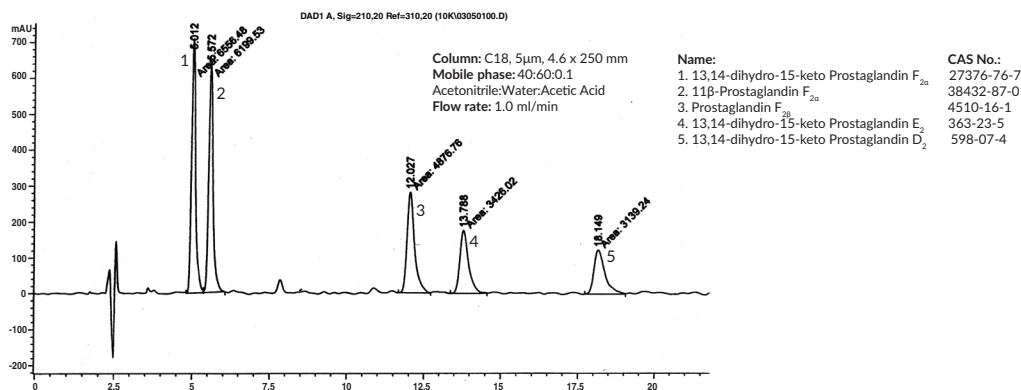
Item No. 10005

Synonym: PG Metabolite HPLC Mixture
Purity: $\geq 98\%$ 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

The prostaglandin metabolite HPLC mixture contains the primary metabolites of five commonly encountered cyclooxygenase products supplied in methyl acetate. This HPLC mixture contains 100 μg of the following prostaglandin metabolites: 13,14-dihydro-15-keto Prostaglandin $\text{F}_{2\alpha}$, 11 β -Prostaglandin $\text{F}_{2\alpha}$, Prostaglandin $\text{F}_{2\beta}$, 13,14-dihydro-15-keto Prostaglandin E_2 , and 13,14-dihydro-15-keto Prostaglandin D_2 .



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, 08/04/2022

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