

# PRODUCT INFORMATION



## Secondary Prostaglandin Metabolite MaxSpec® LC-MS Mixture Item No. 19422

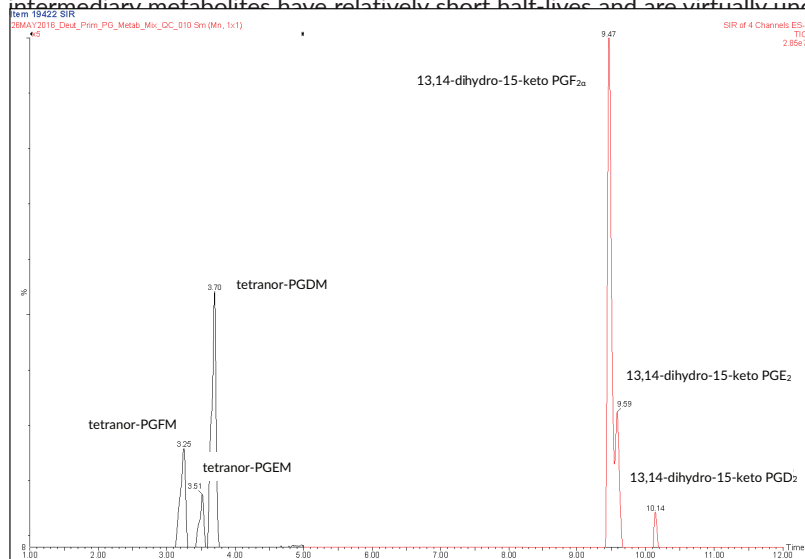
**Purity:** ≥98% for each compound  
**Supplied as:** A solution in acetonitrile (1 µg/ml of each compound)  
**Fill Volume:** >1 ml  
**Storage:** -80°C  
**Stability:** ≥5 years

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

### Description and Contents

This mixture contains the intermediary and secondary metabolites of prostaglandins (PGs) PGE<sub>2</sub> (Item No. 14010), PGD<sub>2</sub> (Item No. 12010), and PGF<sub>2α</sub> (Item No. 16010). The mixture is supplied in an amber ampule in which the headspace has been purged with argon to prevent lipid oxidation. This product has been designed for direct use in LC-MS applications. The solution may be serially diluted for preparation of calibrators and QC standards and/or used directly as a system suitability standard or tuning standard. After opening, we recommend that the mixture be transferred immediately to a 1 ml glass screw cap vial, to prevent solvent evaporation, and stored at -20°C. The mixture should be discarded after multiple freeze/thaw cycles.

The tetranor-PG metabolites represented in this mixture are the major urinary metabolites of PGE<sub>2</sub>, PGD<sub>2</sub>, and PGF<sub>2α</sub> and, while biologically inactive, are used extensively as biomarkers to assess endogenous production of their respective parent PGs. 13,14-dihydro-15-keto PGs are intermediary metabolites of their respective parent PGs formed through the 15-hydroxy PGDH pathway. Though biologically active, these intermediary metabolites have relatively short half-lives and are virtually undetectable in urine.



Item Number: 19422		Secondary Primary Prostaglandin Metabolite MaxSpec® LC-MS Mixture	
Item Number	Item Name	Formula Weight:	MS/MS Transition:
16840	tetranor-PGFM	330.4	329>311
14840	tetranor-PGEM	328.4	327>309
12850	tetranor-PGDM	328.4	327>309
16670	13,14-dihydro-15-keto Prostaglandin F <sub>2α</sub>	354.5	353>193
14650	13,14-dihydro-15-keto Prostaglandin E <sub>2</sub>	352.5	351>113
12610	13,14-dihydro-15-keto Prostaglandin D <sub>2</sub>	352.5	351>175

LC-MS Conditions: Waters Acquity UPLC-Xevo TQ-Smicro  
Mobile Phase A: Water + 0.1% Formic Acid  
Mobile Phase B: Acetonitrile + 0.1% Formic Acid  
Column: Waters BEH C8, 2.1 x 100 mm, 1.7 µm  
Flow Rate: 400 µl/min  
Negative Electrospray Ionization  
SIR Scan

#### 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, 07/14/2023

#### 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