

PRODUCT INFORMATION



Deuterated Primary COX and LOX MaxSpec® LC-MS Mixture Item No. 19228

Supplied as: A solution in ethanol (1 µg/ml of each compound)
Fill volume: >1 ml
Storage: -20°C
Stability: ≥10 years

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

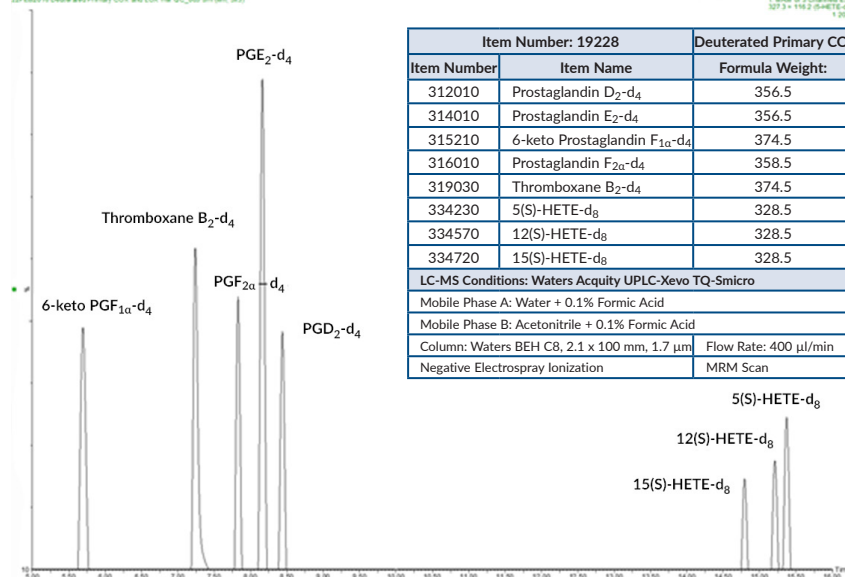
Description

This mixture contains deuterated versions of selected eicosanoids produced by the metabolism of arachidonic acid via the cyclooxygenase and lipoxygenase pathways. 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. This mixture is ideally suited for use as an internal standard in mass spectrometry applications. It should be noted that the mixture contains isobaric analytes. 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 lipid mediators represented in this mixture are produced on demand by various cells and tissue types in response to external stimulation and act locally in an autocrine or paracrine manner. In the body, particularly with thromboxanes and prostaglandins, they are enzymatically converted to inactive metabolites and excreted in the urine. As a result, serum-free media from cultured cells and tissues will serve as the primary sample types for their analysis, although additional sample types could also be used.

Contents

COX and LOX GC MRM
20180215 Deuterated Primary COX and LOX mix GC_MRM (5m, 3s)



1. MRM of 3 Channels (5-
1273 > 116.2 (5-HETE-d8)
1.304

Item Number: 19228		Deuterated Primary COX and LOX MaxSpec® LC-MS Mixture	
Item Number	Item Name	Formula Weight:	MS/MS Transition:
312010	Prostaglandin D ₂ -d ₄	356.5	355>233
314010	Prostaglandin E ₂ -d ₄	356.5	355>319
315210	6-keto Prostaglandin F _{1α} -d ₄	374.5	369>249
316010	Prostaglandin F _{2α} -d ₄	358.5	357>197
319030	Thromboxane B ₂ -d ₄	374.5	373>173
334230	5(S)-HETE-d ₈	328.5	327>116
334570	12(S)-HETE-d ₈	328.5	327>182
334720	15(S)-HETE-d ₈	328.5	327>184

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
MRM 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