

# PRODUCT INFORMATION



## Leukotriene B<sub>4</sub> Pathway MaxSpec<sup>®</sup> LC-MS Mixture Item No. 22640

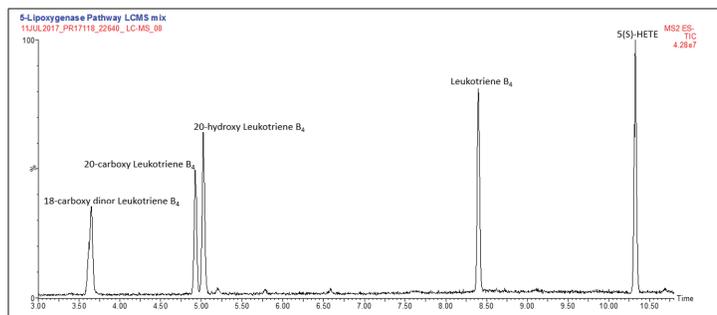
**Supplied as:** A solution in ethanol (1 µg/ml of each compound)  
**Fill Volume:** >1 ml  
**Storage:** -20°C  
**Stability:** ≥5 years

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

### Description and Contents

The Leukotriene B<sub>4</sub> Pathway MaxSpec<sup>®</sup> LC-MS Mixture contains 5-lipoxygenase pathway metabolites derived from arachidonic acid (Item No. 90010). 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 used 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.

This mixture contains 18-carboxy dinor leukotriene B<sub>4</sub> (Item No. 20170), 20-carboxy leukotriene B<sub>4</sub> (Item No. 20180), 20-hydroxy leukotriene B<sub>4</sub> (Item No. 20190), leukotriene B<sub>4</sub> (Item No. 20110), and 5(S)-HETE (Item No. 34230).



Item Number: 22640		Leukotriene B <sub>4</sub> Pathway MaxSpec <sup>®</sup> LC-MS Mixture				
Item Number	Item Name	Formula	Mass (Da)	[MH] <sup>+</sup> (m/z)	RT (min)	
20170	18-carboxy dinor Leukotriene B <sub>4</sub>	C <sub>18</sub> H <sub>26</sub> O <sub>6</sub>	338.4	337.3	3.65	
20180	20-carboxy Leukotriene B <sub>4</sub>	C <sub>20</sub> H <sub>30</sub> O <sub>6</sub>	366.5	365.4	4.92	
20190	20-hydroxy Leukotriene B <sub>4</sub>	C <sub>20</sub> H <sub>32</sub> O <sub>5</sub>	352.5	351.4	5.02	
20110	Leukotriene B <sub>4</sub>	C <sub>20</sub> H <sub>32</sub> O <sub>4</sub>	336.5	335.4	8.40	
34230	5(S)-HETE	C <sub>20</sub> H <sub>32</sub> O <sub>3</sub>	320.5	319.3	10.32	
LC-MS Instrument: Waters Acquity UPLC-Xevo TQD						
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; Injection Volume: 5 µl			
Sample Preparation: 200 µl of the mixture was dried under nitrogen and reconstituted in 200 µl 1:1 ACN:Water						
Negative Electrospray Ionization			Full Mass Scan: 317-369 m/z			

#### 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, 07/17/2023

#### 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