



Certificate of Analysis

Arachidonic Acid Oxylin LC-MS Mixture

Item #: 20666
Batch #: 0652612

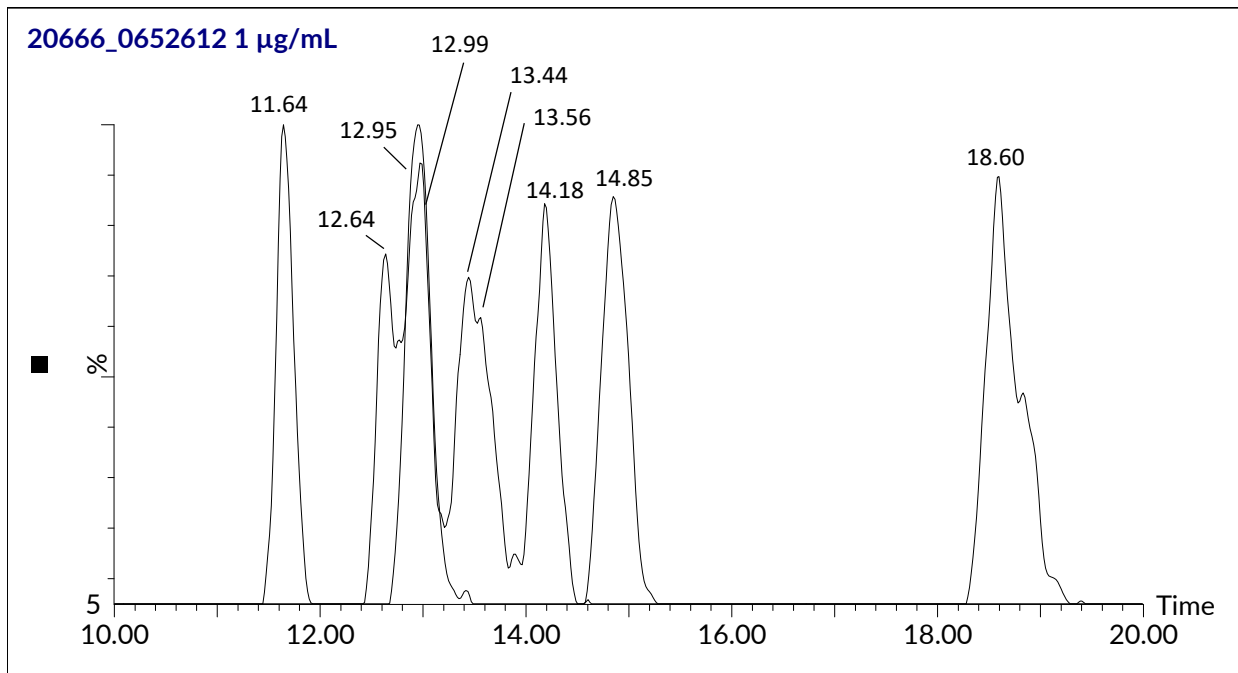
Expiry: 17OCT2027 (valid from date of certification)

Contents: 10 µg/mL (nominal) each in ethanol of 5(S)-HETE, 5-oxoETE, 8(S)-HETE, 9(R)-HETE, 11(S)-HETE, 12(S)-HETE, 12-oxoETE, 15(S)-HETE, 15-oxoETE

Storage and Handling: Temperature sensitive. Product should be stored as supplied at -80°C.

Results:

Method	Specification	Specification Met
Negative ESI LC-MS/MS	All analytes identified	Yes



Cayman Chemical Co.

1180 E. Ellsworth Rd
Ann Arbor, MI 48108
www.caymanchem.com

Contact Us:

1-800-364-9897 or (734)971-3335
Fax: (734)971-3640
Email: techserv@caymanchem.com

For R&D purposes only. Not intended for human or animal use.

Item: 20666 Batch: 0652612		Arachidonic Acid Oxylin LC-MS Mixture		
Item #	Name	Formula Weight	Transitions (m/z)	RT (min)
34720	15(S)-HETE	320.5	319>219	11.64
34510	11(S)-HETE	320.5	319>167	12.64
34730	15-OxoETE	318.5	317>113	12.95
34360	8(S)-HETE	320.5	319>155	12.99
34570	12(S)-HETE	320.5	319>179	13.44
34405	9(R)-HETE	320.5	319>139	13.56
34230	5(S)-HETE	320.5	319>115	14.18
34580	12-OxoETE	318.5	317>153	14.85
34250	5-OxoETE	318.5	317>203	18.60
LC-MS Conditions: Waters Acquity I UPLC - Xevo TQD				
Mobile Phase: A: Water + 0.1% Formic Acid				
Mobile Phase: B: Acetonitrile + 0.1% Formic Acid				
Column: Waters BEH C8, 2.1x100mm, 1.7µm		Flow Rate: 400µL/min		
LC Gradient: Isocratic 48%B over 20 minutes				
Negative Electrospray Ionization				

Mixture was prepared using gravimetric and volumetric solution preparation procedures from a combination of neat and formulated pure standards. The concentration of this mixture was prepared with an estimated error of ±10%.

Cayman Chemical certifies that this product meets the specifications stated in this certificate and warrants this product to meet the stated acceptance criteria through the expiration date when stored unopened as recommended.

Approval:  **Title: Director of Analytical Chemistry**

Certification Date: 17OCT2022