PRODUCT INFORMATION



Lipoxin A₄-d₅ MaxSpec[®] Standard Item No. 24936

CAS Registry No.:	1622429-53-1		
Formal Name:	5S,6R,15S-trihydroxyicosa-7E,9E,11Z,13E	- но	OH
	tetraenoic-19,19,20,20,20-d₅ acid		
Synonyms:	5(S),6(R),15(S)-TriHETE-d ₅ ,		
	$5(S),6(R)-LXA_4-d_5, LXA_4-d_5$		
MF:	$C_{20}H_{27}D_5O_5$		
FW:	357.5		
Purity:	≥95%	ОН	00
Supplied as:	A solution in ethanol; in a deactivated glass ampule		
Concentration:	$10 \mu g/ml$ (nominal); see certificate of analysis for verified concentration		
Storage:	-80°C		
Stability:	≥5 years; Stability testing is ongoing to ensure concentration accuracy. The certificate of analysis and		
	product expiry date will be updated upon completion of testing.		
Special Conditions: Store upright and unopened at -80°C. Warm to room temperature prior to opening			
	Light sensitive.		

Description

Lipoxin A4-d5 (LXA4-d5) is intended for use as an internal standard for the quantification of LXA₄ (Item Nos. 90410 | 10007271) by GC- or LC-mass spectrometry (MS). LXA₄ is a trihydroxy fatty acid containing a conjugated tetraene, produced by the metabolism of 15-HETE or 15-HPETE with human leukocytes.¹ LXA₄ is equipotent to leukotriene B₄ (LTB₄; Item Nos. 20110 | 10007240) in inducing superoxide generation in human neutrophils at 0.1 μ M.² LXA₄ is associated with several other biological functions including leukocyte activation, chemotaxis effects, natural killer cell inhibition, and monocyte migration and adhesion.²⁻⁴

LXA₄-d₅ MaxSpec[®] standard is a quantitative grade standard of LXA₄-d₅ (Item No. 10007737) that has been prepared specifically for mass spectrometry and related applications where quantitative reproducibility is required. The solution has been prepared gravimetrically and is supplied in a deactivated glass ampule sealed under argon. The concentration was verified by comparison to an independently prepared calibration standard. The verified concentration is provided on the certificate of analysis. This LXA_4 -d₅ MaxSpec[®] standard is guaranteed to meet identity, purity, stability, and concentration specifications and is provided with a batch-specific certificate of analysis. Ongoing stability testing is performed to ensure the concentration remains accurate throughout the shelf life of the product. Note: The amount of solution added to the vial is in excess of the listed amount. Therefore, it is necessary to accurately measure volumes for preparation of calibration standards. Follow recommended storage and handling conditions to maintain product quality.

References

- 1. Serhan, C.N., Nicolaou, K.C., Webber, S.E., et al. Lipoxin A. Stereochemistry and biosynthesis. J. Biol. Chem. 261(35), 16340-16345 (1986).
- 2 Serhan, C.N., Hamberg, M., and Samuelsson, B. Lipoxins: Novel series of biologically active compounds formed from arachidonic acid in human leukocytes. Proc. Natl. Acad. Sci. USA 81(17), 5335-5339 (1984).
- 3 Ramstedt, U., Serhan, C.N., Nicolaou, K.C., et al. Lipoxin A-induced inhibition of human natural killer cell cytotoxicity: Studies on stereospecificity of inhibition and mode of action. J. Immunol. 138(1), 266-270 (1987).
- 4. Maddox, J.F., and Serhan, C.N. Lipoxin A_{4} and B_{4} are potent stimuli for human monocyte migration and adhesion: Selective inactivation by dehydrogenation and reduction. J. Exp. Med. 183(1), 137-146 (1996).

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

uyer 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, 06/21/2021

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