

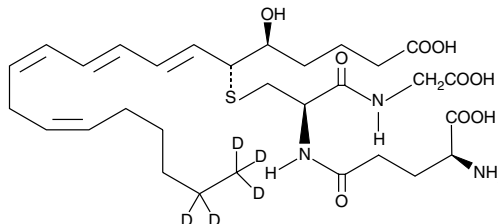
Product Information



Leukotriene C₄-d₅

Item No. 10006198

CAS Registry No.: 1441421-73-3
Formal Name: 5S-hydroxy-6R-(S-glutathionyl)-7E,9E,11Z,14Z-d₅-eicosatetraenoic acid
Synonym: LTC₄-d₅
MF: C₃₀H₄₂D₅N₃O₉S
FW: 630.8
Purity: ≥97%
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
Stability: ≥1 year at -80°C
Supplied as: A solution in ethanol:water (95:5)
UV/Vis.: λ_{max}: 280 nm ε: 40,000
Miscellaneous: Light Sensitive



Laboratory Procedures

Leukotriene C₄-d₅ (LTC₄-d₅) contains five deuterium atoms at the 19, 19', 20, 20 and 20 positions. It is intended for use as an internal standard for the quantification of LTC₄ by GC- or LC-mass spectrometry (MS). For long term storage, we suggest that LTC₄-d₅ be stored as supplied at -80°C. It will be stable for at least one year.

LTC₄-d₅ is supplied as a solution in ethanol:water (95:5). To change the solvent, simply evaporate the ethanol:water under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of LTC₄-d₅ in these solvents is approximately 50 mg/ml.

LTC₄-d₅ is used as an internal standard for the quantification of LTC₄ by stable isotope dilution MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated *versus* unlabeled).

LTC₄ is the parent cysteinyl leukotriene produced by the addition of glutathione to LTA₄. LTC₄ is produced by neutrophils, macrophages, and mast cells, and by *trans*-cellular metabolism in platelets.¹ It is one of the constituents of slow-reacting substance of anaphylaxis (SRS-A) and exhibits potent smooth muscle contracting activity.² LTC₄-induced bronchoconstriction and enhanced vascular permeability contribute to the pathogenesis of asthma and acute allergic hypersensitivity.^{3,4} The concentration of LTC₄ required to produce marked contractions of lung parenchymal strips and isolated tracheal rings is about 1 nM.⁴

References

1. Maclouf, J.A. and Murphy, R.C. Transcellular metabolism of neutrophil-derived leukotriene A₄ by human platelets. A potential cellular source of leukotriene C₄. *J. Biol. Chem.* **263**, 174-181 (1988).
2. Piper, P.J. Formation and actions of leukotrienes. *Physiol. Rev.* **64**, 744-761 (1984).
3. Samuelsson, B. Leukotrienes: Mediators of immediate hypersensitivity reactions and inflammation. *Science* **220**, 568-575 (1983).
4. Lefer, A.M. Leukotrienes as mediators of ischemia and shock. *Biochem. Pharmacol.* **35**, 123-127 (1986).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10006198

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY; NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog.**

Copyright Cayman Chemical Company, 12/12/2013

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com