

# PRODUCT INFORMATION



## Leukotriene E<sub>4</sub>-d<sub>5</sub> MaxSpec<sup>®</sup> Standard

Item No. 25299

CAS Registry No.: 1240398-14-4

Formal Name: 5S-hydroxy-6R-(S-cysteinyl)-  
7E,9E,11Z,14Z-eicosatetraenoic-  
19,19,20,20,20-d<sub>5</sub> acid

Synonym: LTE<sub>4</sub>-d<sub>5</sub>

MF: C<sub>23</sub>H<sub>32</sub>D<sub>5</sub>NO<sub>5</sub>S

FW: 444.6

Purity: ≥95%

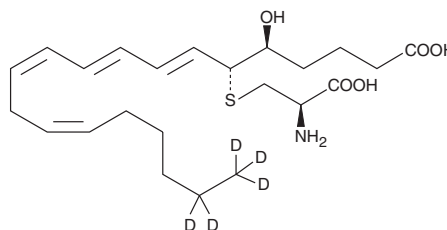
Supplied as: A solution in ethanol; in a deactivated glass ampule

Concentration: 10 µg/ml (nominal); see certificate of analysis for verified concentration

Storage: -80°C

Stability: ≥7 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

Leukotriene E<sub>4</sub>-d<sub>5</sub> (LTE<sub>4</sub>-d<sub>5</sub>) is intended for use as an internal standard for the quantification of LTE<sub>4</sub> by GC- or LC-mass spectrometry. LTE<sub>4</sub> is produced by the action of dipeptidase on LTD<sub>4</sub>, leaving only the cysteinyl group still attached to the fatty acid backbone.<sup>1</sup> It is one of the constituents of slow-reacting substance of anaphylaxis (SRS-A).<sup>2</sup> LTE<sub>4</sub> is considerably less active (8 to 12-fold) than LTC<sub>4</sub> in the biological activities characteristic of cysteinyl leukotrienes.<sup>1,3</sup> Unlike LTC<sub>4</sub> and LTD<sub>4</sub>, LTE<sub>4</sub> accumulates in both plasma and urine. Therefore, urinary excretion of LTE<sub>4</sub> is most often used as an indicator of asthma.<sup>4-6</sup> In humans, basal levels of LTE<sub>4</sub> range from 1-100 pg/mg creatinine. In asthmatic patients, urinary LTE<sub>4</sub> levels increase to 80-1,000 pg/mg creatinine.<sup>5</sup>

LTE<sub>4</sub>-d<sub>5</sub> MaxSpec<sup>®</sup> standard is a quantitative grade standard of LTE<sub>4</sub>-d<sub>5</sub> (Item No. 10007858) that has been prepared specifically for mass spectrometry or any application 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. This LTE<sub>4</sub>-d<sub>5</sub> MaxSpec<sup>®</sup> 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. Bernström, K. and Hammarström, S. *J. Biol. Chem.* **256**(18), 9579-9582 (1981).
2. Samuelsson, B. *Science* **220**(4597), 568-575 (1983).
3. Lefer, A.M. *Biochem. Pharmacol.* **35**(2), 123-127 (1986).
4. Kumlin, M., Stensvad, F., Larsson, L., et al. *Clin. Exp. Allergy* **25**(5), 467-479 (1995).
5. Drazen, J.M., O'Brien, J., Sparrow, D., et al. *Am. Rev. Respir. Dis.* **146**(1), 104-108 (1992).
6. Kumlin, M., Dahlén, B., Björck, T., et al. *Am. Rev. Respir. Dis.* **146**(1), 96-103 (1992).

#### 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, 01/24/2023

#### 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