

PRODUCT INFORMATION



MCTR1

Item No. 17007

CAS Registry No.: 1784701-61-6
Formal Name: L-γ-glutamyl-S-[(1R,2E,4E,6Z,9Z)-12-carboxy-1-[(1S,3Z,6Z)-1-hydroxy-3,6-nonadien-1-yl]-2,4,6,9-dodecatetraen-1-yl]-L-cysteinyl-glycine
Synonyms: 13-glutathionyl-14-hydroxy Docosahexaenoic Acid, Maresin Conjugates in Tissue Regeneration 1, Maresin Sulfido Conjugate 1

MF: C₃₂H₄₇N₃O₉S

FW: 649.8

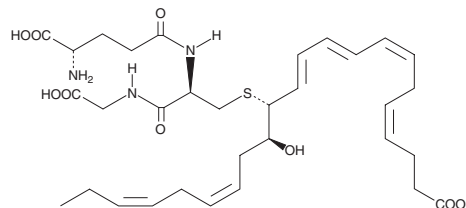
Purity: ≥95%

UV/Vis.: λ_{max}: 282 nm

Supplied as: A solution in ethanol

Storage: -80°C

Stability: ≥1 year



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

Laboratory Procedures

Maresin conjugates in tissue regeneration 1 (MCTR1) is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of MCTR1 in ethanol is approximately 1 mg/ml and approximately 50 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. If an organic solvent-free solution of MCTR1 is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of MCTR1 in PBS, pH 7.2, is approximately 0.1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

MCTR1 is a specialized pro-resolving mediator (SPM) synthesized from docosahexaenoic acid (DHA; Item No. 90310) in macrophages at the site of inflammation.^{1,2} DHA is oxidized to maresin 1 (MaR1; Item No. 10878), which is then converted to MCTR1 by glutathione S-transferase Mu 4 or leukotriene C₄ synthase.³⁻⁵ MCTR1 accelerates tissue regeneration in planaria (1 and 100 nM).² Pretreatment with MCTR1 (50 ng/mouse, i.p.) prior to *E. coli* administration reduces neutrophil infiltration, shortens the inflammatory resolution period, and increases phagocytosis of *E. coli* by macrophages.² When administered at a dose of 100 ng 12h post *E. coli* infection in a mouse model of peritonitis, MCTR1 reduces the amount of eicosanoids in the exudate.

References

1. Serhan, C.N. *Nature* **510(7503)**, 92-101 (2014).
2. Dalli, K.J., Sanger, J.M., Rodriguez, A.R., et al. *PLoS One* **11(2)**, e0149319 (2016).
3. Dalli, J., Chiang, N., and Serhan, C.N. *Proc. Natl. Acad. Sci. USA* **111(44)**, E4753-E4761 (2014).
4. Dalli, J., Vlasakov, I., Riley, I.R., et al. *Proc. Natl. Acad. Sci. USA* **113(43)**, 12232-12237 (2016).
5. Serhan, C.N., Dalli, J., Karamnov, S., et al. *FASEB J.* **26(4)**, 1755-1765 (2012).

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, 10/31/2022

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