

PRODUCT INFORMATION



Resolvin D3-d₅ Item No. 19512

Formal Name: 4S,11R,17S-trihydroxydocosa-5Z,7E,9E,13Z,15E,19Z-hexaenoic-21,21,22,22,22-d₅ acid

Synonym: RVD3-d₅

MF: C₂₂H₂₇D₅O₅

FW: 381.5

Chemical Purity: ≥98% (Resolvin D3)

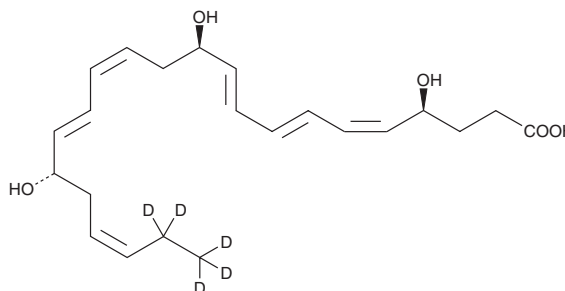
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀

UV/Vis.: λ_{max}: 240, 262, 273, 283 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

Resolvin D3-d₅ (RVD3-d₅) is intended for use as an internal standard for the quantification of RVD3 (Item No. 13834) by GC- or LC-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).

RVD3-d₅ 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. It is recommended that this product be stored and handled in an ethanol solution. Resolvins can isomerize and degrade when put into freeze thaw conditions and/or in solvents such as dimethyl formamide or DMSO.

Description

RVD3 is a DHA-derived product first identified in mouse inflammatory exudates.¹ It reduces neutrophil infiltration *in vivo* in both mouse peritonitis and dermal inflammation. In addition to suppressing leukocyte migration, RVD3 enhances macrophage phagocytosis and efferocytosis.² Unlike other resolvins, RVD3 appears late in resolution in mouse peritonitis.^{2,3}

References

1. Serhan, C.N., Hong, S., Gronert, K., *et al.* *J. Exp. Med.* **196**(8),1025-1037 (2002).
2. Dalli, J., Winkler, J.W., Colas, R.A., *et al.* *Chem. Biol.* **20**(2), 188-201 (2015).
3. Serhan, C.N. *Nature* **510**(7503), 92-101 (2014).

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, 06/18/2019

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