

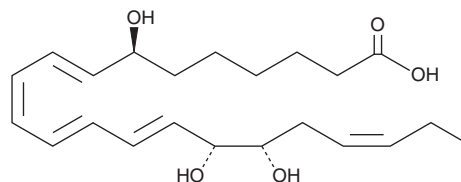
PRODUCT INFORMATION



Resolvin D2 n-3 DPA

Item No. 34482

CAS Registry No.: 2093111-29-4
Formal Name: 7S,16R,17S-trihydroxy-8E,10Z,12E,14E,19Z-docosapentaenoic acid
Synonyms: 7(S),16(R),17(S)-Resolvin D2 n-3 DPA, RvD2 n-3 DPA
MF: C₂₂H₃₄O₅
FW: 378.5
Purity: ≥98%
UV/Vis.: λ_{max}: 289, 302, 317 nm
Supplied as: A 50 µg/ml 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.

Description

Resolvin D2 n-3 DPA (RvD2 n-3 DPA) is a specialized pro-resolving mediator (SPM).¹ It is formed from docosapentaenoic acid (n-3 DPA; Item Nos. 90165 | 21907), an intermediate in the conversion of eicosapentaenoic acid (Item Nos. 90110 | 90110.1 | 21908) to docosahexaenoic acid (Item Nos. 90310 | 17950), in human leukocytes. RvD2 n-3 DPA (1 nM) reduces TNF-α-induced chemotaxis and adhesion of isolated human neutrophils. *In vivo*, RvD2 n-3 DPA (100 ng/animal; i.v.) reduces peritoneal neutrophil infiltration and exudate levels of IL-6 and chemokine (C-C motif) ligand 2 (CCL2) in a mouse model of zymosan-induced inflammation.

Reference

1. Dalli, J., Colas, R.A., and Serhan, C.N. Novel n-3 immunoresolvents: Structures and actions. *Sci. Rep.* **3**, 1940 (2013).

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, 01/14/2026

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