

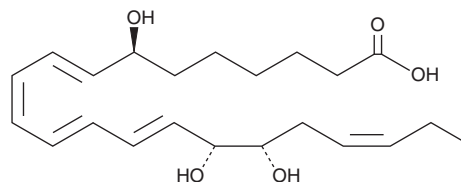
# PRODUCT INFORMATION



## 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:** C<sub>22</sub>H<sub>34</sub>O<sub>5</sub>  
**FW:** 378.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 289, 302, 317 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.

### Description

Resolvin D2 n-3 DPA (RvD2 n-3 DPA) is a specialized pro-resolving mediator (SPM).<sup>1</sup> 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, 08/19/2021

#### 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