

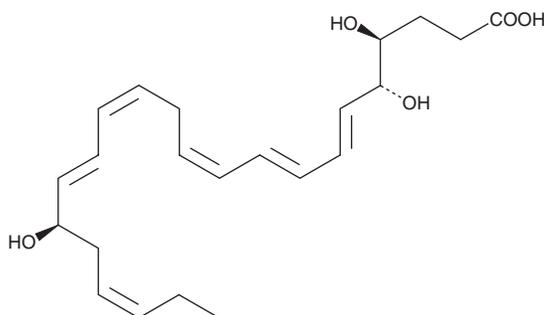
# PRODUCT INFORMATION



## 17(R)-Resolvin D4

Item No. 9002881

**CAS Registry No.:** 528583-89-3  
**Formal Name:** 4S,5R,17R-trihydroxy-6E,8E,10Z,13Z,15E,19Z-docosahexaenoic acid  
**Synonyms:** 4(S),5(R),17(R)-trihydroxy-DHA, 17(R)-RvD4  
**MF:** C<sub>22</sub>H<sub>32</sub>O<sub>5</sub>  
**FW:** 376.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 228, 277, 332 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

17(R)-Resolvin D4 is supplied as a solution in ethanol. To change the solvent, simply evaporate the 17(R)-RvD4 under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol and dimethyl formamide purged with an inert gas can be used. The solubility of 17(R)-RvD4 in these solvents is approximately 50 mg/ml.

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 17(R)-RvD4 is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of 17(R)-resolvin D4 in PBS, pH 7.2, is approximately 0.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

17(R)-RvD4 is an aspirin-triggered epimer of RvD4 (Item No. 13835).<sup>1</sup>

### Reference

1. Serhan, C.N., Hong, S., Gronert, K., *et al.* Resolvins: A family of bioactive products of ω-3 fatty acid transformation circuits by aspirin treatment that counter proinflammation signals. *J. Exp. Med.* **196(8)**, 1025-1037 (2002).

**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, 11/30/2017

### 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