

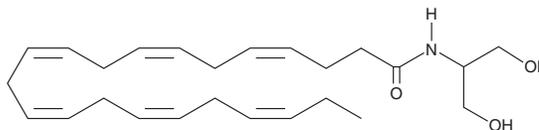
# PRODUCT INFORMATION



## Docosahexaenoyl Serinol

Item No. 35678

**CAS Registry No.:** 1638355-62-0  
**Formal Name:** N-(1,3-dihydroxypropan-2-yl) docosa-4Z,7Z,10Z,13Z,16Z,19Z-hexaenamide  
**Synonym:** Docosahexaenoyl Serinolamide  
**MF:** C<sub>25</sub>H<sub>39</sub>NO<sub>3</sub>  
**FW:** 401.6  
**Purity:** ≥95%  
**Supplied as:** A 5 mg/ml solution in ethanol  
**Storage:** 20°C  
**Stability:** ≥1 year



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

### Laboratory Procedures

Docosahexaenoyl serinol 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. Docosahexaenoyl serinol is soluble in chloroform.

### Description

Docosahexaenoyl serinol is a derivative of 2-docosahexaenoyl glycerol (2-DG). It inhibits the growth of A549 human lung adenocarcinoma cells (IC<sub>50</sub> = 0.48 μM).<sup>1</sup> It also inhibits degranulation of RBL-2H3 mast cells (IC<sub>50</sub> = 16 μM).<sup>2</sup> Oral administration of docosahexaenoyl serinol (20 mg/animal) suppresses the IgE-dependent passive cutaneous anaphylaxis (PCA) reaction by 17% in mice.

### References

1. Tremblay, H., St-Georges, C., Legault, M.A., *et al.* One-pot synthesis of polyunsaturated fatty acid amides with anti-proliferative properties. *Bioorg. Med. Chem. Lett.* **24(24)**, 5635-5638 (2014).
2. Kim, I.-H., Kanayama, Y., Nishiwaki, H., *et al.* Structure-activity relationships of fish oil derivatives with anti-allergic activity in vitro and in vivo. *J. Med. Chem.* **62(21)**, 9576-9592 (2019).

#### 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, 01/26/2026

#### 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