

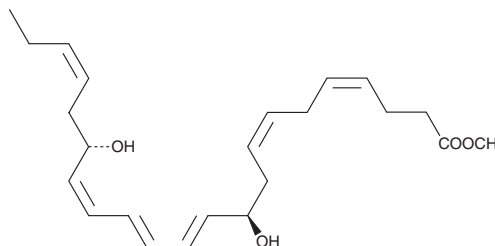
# PRODUCT INFORMATION



## Protectin D1 methyl ester

Item No. 36746

**CAS Registry No.:** 934005-36-4  
**Formal Name:** 10R,17S-dihydroxy-4Z,7Z,11E,13E,15Z,19Z-docosahexaenoic acid, methyl ester  
**Synonyms:** Neuroprotectin D1 methyl ester, NPD1 methyl ester, PD1 methyl ester  
**MF:** C<sub>23</sub>H<sub>34</sub>O<sub>4</sub>  
**FW:** 374.5  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 271, 282 nm  
**Supplied as:** A 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.

### Description

Protectin D1 methyl ester is an esterified form of the specialized pro-resolving mediator protectin D1 (Item No. 10010390). It is an intermediate in the synthesis of protectin D1.<sup>1</sup> Protectin D1 methyl ester (333 µg/kg) reduces infarct area and volume and increases blood vessel density in the tissue surrounding the infarct in a rat model of ischemic stroke induced by middle cerebral artery occlusion (MCAO).<sup>2</sup>

### References

1. Rodriguez, A.R. and Spur, B.W. Total synthesis of the potent anti-inflammatory lipid mediator Protectin D1. *Tetrahedron Lett.* **55**(43), 6011-6015 (2014).
2. Bazan, N.G., Eady, T.N., Khoutorova, L., *et al.* Novel aspirin-triggered neuroprotectin D1 attenuates cerebral ischemic injury after experimental stroke. *Exp. Neurol.* **236**(1), 122-130 (2012).

#### 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, 04/25/2022

#### 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](http://WWW.CAYMANCHEM.COM)