

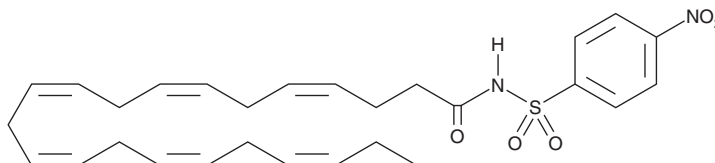
PRODUCT INFORMATION



pNPS-DHA

Item No. 29507

CAS Registry No.: 2454246-25-2
Formal Name: (4Z,7Z,10Z,13Z,16Z,19Z)-N-[(4-nitrophenyl)sulfonyl]docosa-4,7,10,13,16,19-docosaheptaenamide
Synonym: CAY10741, Compound 59, DHA-nitrophenyl sulfonamide, p-Nitrophenyl sulfonamide, Docosaheptaenoic Acid
MF: C₂₈H₃₆N₂O₅S
FW: 512.7
Purity: ≥98%
UV/Vis.: λ_{max}: 256 nm
Supplied as: A 10 mg/ml solution in methyl acetate
Storage: -20°C
Stability: ≥4 years



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

Description

pNPS-DHA is an arylamide derivative of docosaheptaenoyl ethanolamide (DHEA; Item No. 10007534) that has anti-allergic activity.¹ It inhibits degranulation of RBL-2H3 mast cells (IC₅₀ = 15 μM). pNPS-DHA (1,000 mg/kg) inhibits IgE-dependent passive cutaneous anaphylaxis (PCA) in mice.

Reference

1. 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, 07/10/2026

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