

PRODUCT INFORMATION



H2L5186303

Item No. 14663

CAS Registry No.: 139262-76-3

Formal Name: (Z,Z)-4,4'-[1,3-phenylenebis(oxy-4,1-phenyleneimino)]bis[4-oxo-2-butenic acid]

MF: C₂₆H₂₀N₂O₈

FW: 488.5

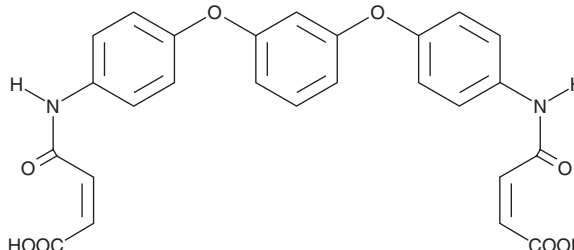
Purity: ≥98%

UV/Vis.: λ_{max}: 298 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

H2L5186303 is supplied as a crystalline solid. A stock solution may be made by dissolving the H2L5186303 in the solvent of choice, which should be purged with an inert gas. H2L5186303 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of H2L5186303 in these solvents is approximately 15 and 20 mg/ml, respectively.

H2L5186303 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, H2L5186303 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. H2L5186303 has a solubility of approximately 0.11 mg/ml in a 1:8 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

H2L5186303 is a selective lysophosphatidic acid 2 (LPA₂) receptor antagonist (IC₅₀ = 9 nM in a LPA-elicited calcium mobilization assay).¹ It inhibits LPA₁ and LPA₃ at much higher concentrations (IC₅₀s = 27,354 and 4,504 nM, respectively).¹

Reference

1. Fells, J. I., Tsukahara, R., Liu, J. *et al.* Structure-based drug design identifies novel LPA₃ antagonists. *Bioorg. Med. Chem.* **17**(21), 7457-7464 (2009).

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, 11/01/2022

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