

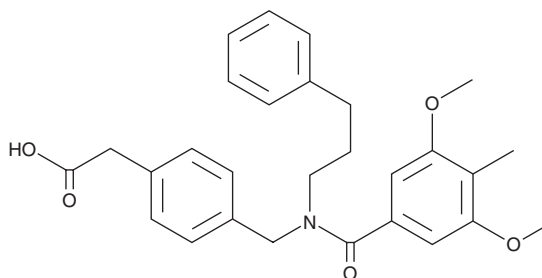
PRODUCT INFORMATION

ONO-7300243

Item No. 22052



CAS Registry No.: 638132-34-0
Formal Name: 4-[[[(3,5-dimethoxy-4-methylbenzoyl)(3-phenylpropyl)amino]methyl]-benzeneacetic acid
MF: C₂₈H₃₁NO₅
FW: 461.6
Purity: ≥98%
Supplied as: A 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

ONO-7300243 is supplied as a solid. A stock solution may be made by dissolving the ONO-7300243 in the solvent of choice, which should be purged with an inert gas. ONO-7300243 is soluble in the organic solvent DMSO at a concentration of approximately 50 mg/ml.

Description

ONO-7300243 is an antagonist of lysophosphatidic acid receptor 1 (LPA₁; IC₅₀ = 0.16 μM).¹ It is selective for LPA₁ over the LPA₂ and LPA₃ receptors (IC₅₀s = 8.6 and >10 μM, respectively). ONO-7300243 dose-dependently reduces intraurethral pressure induced by LPA in rats (ID₅₀ = 11.6 mg/kg) without affecting mean blood pressure.

Reference

1. Terkado, M., Suzuki, H., Hashimura, K., *et al.* Discovery of ONO-7300243 from a novel class of lysophosphatidic acid receptor 1 antagonists: From hit to lead. *ACS Med. Chem. Lett.* **7**(10), 913-918 (2016).

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, 12/05/2022

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