

PRODUCT INFORMATION



Prostaglandin D₂ Inhibitor

Item No. 42095

CAS Registry No.: 885066-67-1

Formal Name: 3-[6-[[2-(2,4-dichlorophenyl)ethyl]amino]-2-methoxy-4-pyrimidinyl]- α,α -dimethyl-benzeneacetic acid

Synonyms: PGD₂-IN-1, PGD₂ Inhibitor, PGD₂ Inhibitor 1, Prostaglandin D₂-IN-1, Prostaglandin D₂ Inhibitor 1

MF: C₂₃H₂₃Cl₂N₃O₃

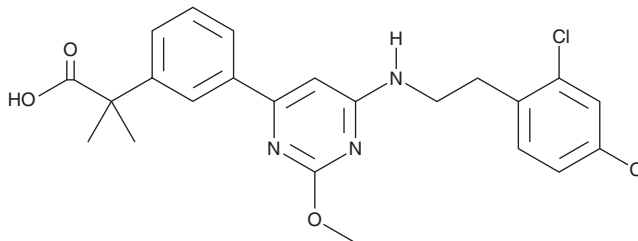
FW: 460.4

Purity: $\geq 98\%$

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

Prostaglandin D₂ (PGD₂) inhibitor is supplied as a solid. A stock solution may be made by dissolving the PGD₂ inhibitor in the solvent of choice, which should be purged with an inert gas. PGD₂ inhibitor is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

PGD₂ inhibitor is an antagonist of PGD₂ receptor (DP; IC₅₀ = 0.3 nM).¹ It inhibits PGD₂-induced cAMP accumulation in LS 174T cells that endogenously express the DP receptor.

Reference

1. Mu, L., Aguiar, J., Ardati, A., *et al.* Understanding DP receptor antagonism using a CoMSIA approach. *Bioorg. Med. Chem. Lett.* **21**(1), 66-75 (2011).

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, 10/31/2024

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