

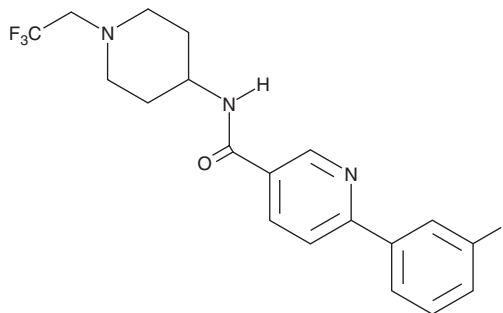
PRODUCT INFORMATION



Prostaglandin D Synthase (hematopoietic-type) Inhibitor I

Item No. 16256

CAS Registry No.: 1033836-12-2
Formal Name: 6-(3-fluorophenyl)-N-[1-(2,2,2-trifluoroethyl)-4-piperidinyl]-3-pyridinecarboxamide
Synonym: H-PGDS Inhibitor I
MF: C₁₉H₁₉F₄N₃O
FW: 381.4
Purity: ≥98%
UV/Vis.: λ_{max}: 260, 287 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

Prostaglandin D synthase (hematopoietic-type) (H-PGDS) inhibitor I is supplied as a crystalline solid. A stock solution may be made by dissolving the H-PGDS inhibitor I in the solvent of choice, which should be purged with an inert gas. H-PGDS inhibitor I is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of H-PGDS inhibitor I in ethanol is approximately 1 mg/ml and approximately 5 mg/ml in DMSO and DMF.

Description

Prostaglandin D synthase (hematopoietic-type) (H-PGDS) inhibitor I is an inhibitor of H-PGDS (IC₅₀ = 32 nM for the human enzyme).¹ It is selective for H-PGDS over lipocalin-type PGDS, microsomal prostaglandin E synthase-1 (mPGES-1), COX-1, COX-2, and 5-lipoxygenase (5-LO; IC₅₀s = >10,000 nM for all). H-PGDS inhibitor I (1 or 10 mg/kg) inhibits PGD₂ (Item No. 12010) production *in vivo* in rat spleen. It completely inhibits allergen-induced airway hyperresponsiveness in an ovine model of asthma.

Reference

1. Carron, C.P., Trujillo, J.I., Olson, K.L., *et al.* Discovery of an oral potent selective inhibitor of hematopoietic prostaglandin D synthase (HPGDS). *ACS Med. Chem. Lett.* **1**(2), 59-63 (2010).

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, 12/08/2022

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