

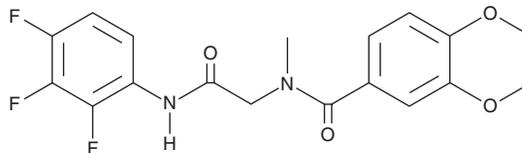
PRODUCT INFORMATION



IMB-808

Item No. 42448

CAS Registry No.: 870768-70-0
Formal Name: 2,3-dihydro-N-methyl-N-[2-oxo-2-[(2,3,4-trifluorophenyl)amino]ethyl]-1,4-benzodioxin-6-carboxamide
MF: C₁₈H₁₅F₃N₂O₄
FW: 380.3
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

IMB-808 is supplied as a solid. A stock solution may be made by dissolving the IMB-808 in the solvent of choice, which should be purged with an inert gas. IMB-808 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

IMB-808 is slightly soluble (0.1-1 mg/ml) in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

IMB-808 is an agonist of liver X receptor α (LXR α) and LXR β .¹ It induces transactivation of LXR α and LXR β in reporter assays using HEK293T cells (EC₅₀s = 0.15 and 0.53 μ M, respectively). IMB-808 (1 and 10 μ M) increases protein levels of ATP-binding cassette transporter 1, subfamily G (ABCG1) and ATP-binding cassette transporter 1, subfamily A (ABCA1) in RAW 264.7 and THP-1 cells. It induces cholesterol efflux and reduces oxidized LDL-induced lipid accumulation in RAW 264.7 cells when used at concentrations of 1 or 10 μ M.

Reference

1. Li, N., Wang, X., Xu, Y., *et al.* Identification of a novel liver X receptor agonist that regulates the expression of key cholesterol homeostasis genes with distinct pharmacological characteristics. *Mol. Pharmacol.* **91**(4), 264-276 (2017).

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, 11/08/2024

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