

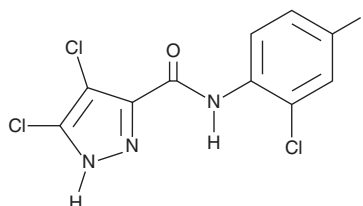
PRODUCT INFORMATION



BLX3887

Item No. 27391

CAS Registry No.: 934758-70-0
Formal Name: 4,5-dichloro-N-(2-chloro-4-fluorophenyl)-1H-pyrazole-3-carboxamide
Synonym: 15-LO-1 Inhibitor 3887
MF: C₁₀H₅Cl₃FN₃O
FW: 308.5
Purity: ≥95%
UV/Vis.: λ_{max}: 260 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

BLX3887 is supplied as a crystalline solid. A stock solution may be made by dissolving the BLX3887 in the solvent of choice, which should be purged with an inert gas. BLX3887 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of BLX3887 in these solvents is approximately 15 and 30 mg/ml, respectively.

BLX3887 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BLX3887 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. BLX3887 has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

BLX3887 is an inhibitor of 15-lipoxygenase type 1 (15-LO-1; IC₅₀ = 32 nM in a cell-free enzyme assay).¹ It is selective for 15-LO-1 over 15-LO-2, which it does not inhibit, 5-LO (IC₅₀ = 472 nM), and 12-LO (IC₅₀ = 3,310 nM). BLX3887 inhibits the production of 15-LO metabolites selectively in eosinophils over neutrophils when used at a concentration of 10 μM. It also inhibits endocytosis in, and the migration of, isolated human peripheral blood mononuclear cell-derived dendritic cells *in vitro*.

Reference

1. Archambault, A.-S., Turcotte, C., Martin, C., *et al.* Comparison of eight 15-lipoxygenase (LO) inhibitors on the biosynthesis of 15-LO metabolites by human neutrophils and eosinophils. *PLoS One* **13**(8), e0202424 (2018).

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, 07/02/2025

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