

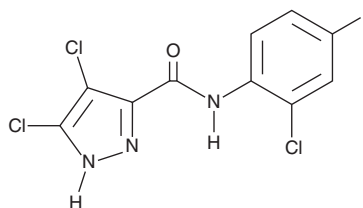
# PRODUCT INFORMATION



**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:** C<sub>10</sub>H<sub>5</sub>Cl<sub>3</sub>FN<sub>3</sub>O  
**FW:** 308.5  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 260 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 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; IC<sub>50</sub> = 32 nM in a cell-free enzyme assay).<sup>1</sup> It is selective for 15-LO-1 over 15-LO-2, which it does not inhibit, 5-LO (IC<sub>50</sub> = 472 nM), and 12-LO (IC<sub>50</sub> = 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, 01/22/2020

## 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