

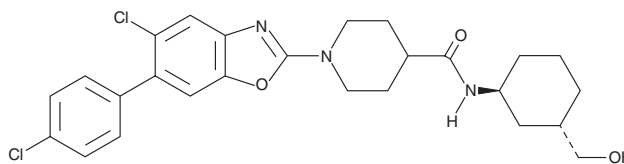
# PRODUCT INFORMATION



**PF-4693627**

Item No. 22768

**CAS Registry No.:** 1312815-93-2  
**Formal Name:** 1-[5-chloro-6-(4-chlorophenyl)-2-benzoxazolyl]-N-[(1S,3S)-3-(hydroxymethyl)cyclohexyl]-4-piperidinecarboxamide  
**MF:** C<sub>26</sub>H<sub>29</sub>Cl<sub>2</sub>N<sub>3</sub>O<sub>3</sub>  
**FW:** 502.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 300 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

PF-4693627 is supplied as a crystalline solid. A stock solution may be made by dissolving the PF-4693627 in the solvent of choice, which should be purged with an inert gas. PF-4693627 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of PF-4693627 in ethanol is approximately 5 mg/ml and approximately 10 mg/ml in DMSO and DMF.

PF-4693627 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PF-4693627 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. PF-4693627 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

PF-4693627 is an orally bioavailable inhibitor of microsomal prostaglandin E (PGE) synthase-1 (mPGES-1; IC<sub>50</sub> = 3 nM).<sup>1</sup> It is selective for mPGES-1 over PGDS, TXAS, and 5-LO in HWB-1483 cells, which have IC<sub>50</sub> values greater than 50 μM, as well as COX-2 in fetal fibroblasts (IC<sub>50</sub> = >10 μM). PF-4693627 inhibits LPS-stimulated synthesis of PGE<sub>2</sub> in human whole blood (IC<sub>50</sub> = 109 nM). *In vivo*, PF-4693627 (10 mg/kg) inhibits PGE<sub>2</sub> production in a guinea pig model of carrageenan-induced air pouch inflammation.

## Reference

1. Arhancet, G.B., Walker, D.P., Metz, S., *et al.* Discovery and SAR of PF-4693627, a potent, selective and orally bioavailable mPGES-1 inhibitor for the potential treatment of inflammation. *Bioorg. Med. Chem. Lett.* 23(4), 1114-1119 (2013).

### 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, 12/19/2022

## 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