

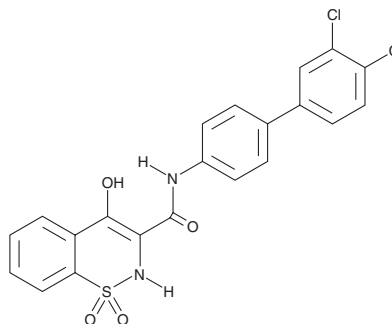
# PRODUCT INFORMATION



## PF-9184

Item No. 18904

**CAS Registry No.:** 1221971-47-6  
**Formal Name:** N-(3',4'-dichloro[1,1'-biphenyl]-4-yl)-4-hydroxy-2H-1,2-benzothiazine-3-carboxamide 1,1-dioxide  
**MF:** C<sub>21</sub>H<sub>14</sub>Cl<sub>2</sub>N<sub>2</sub>O<sub>4</sub>S  
**FW:** 461.3  
**Purity:** ≥98%  
**Supplied as:** A crystalline solid  
**UV/Vis.:** λ<sub>max</sub>: 269, 335 nm  
**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-9184 is supplied as a crystalline solid. A stock solution may be made by dissolving the PF-9184 in the solvent of choice. PF-9184 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of PF-9184 in these solvents is approximately 15 and 20 mg/ml, respectively.

PF-9184 is sparingly soluble 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

Microsomal prostaglandin E synthase-1 (mPGES-1) converts the COX product prostaglandin H<sub>2</sub> (PGH<sub>2</sub>; Item No. 17020) into the biologically active PGE<sub>2</sub> (Item No. 14010).<sup>1</sup> Like COX-2, the expression of mPGES-1 is induced in response to pro-inflammatory mediators, including LPS, IL-1β, and TNF-α.<sup>2</sup> PF-9184 is a potent inhibitor of mPGES-1 (IC<sub>50</sub> = 16.5 nM for recombinant human enzyme).<sup>3</sup> It displays at least 6,500-fold selectivity for mPGES-1 over COX-1 and COX-2. PF-9184 blocks the synthesis of PGE<sub>2</sub> in LPS-treated human whole blood and in IL-1β-stimulated fibroblasts (IC<sub>50</sub>s = 0.4-5 μM).<sup>4,5</sup>

## References

1. Jakobsson, P.-J., Thorén, S., Morgenstern, R., *et al.* Identification of human prostaglandin E synthase: A microsomal, glutathione-dependent, inducible enzyme, constituting a potential novel drug target. *Proc. Natl. Acad. Sci. USA* **96**, 7220-7225 (1999).
2. Stichtenoth, D.O., Thorén, S., Bian, H., *et al.* Microsomal prostaglandin E synthase is regulated by proinflammatory cytokines and glucocorticoids in primary rheumatoid synovial cells. *J. Immunol.* **167**, 469-474 (2001).
3. Mbalaviele, G., Pauley, A.M., Shaffer, A.F., *et al.* Distinction of microsomal prostaglandin E synthase-1 (mPGES-1) inhibition from cyclooxygenase-2 inhibition in cells using a novel, selective mPGES-1 inhibitor. *Biochem. Pharmacol.* **79(10)**, 1445-1454 (2010).
4. Chang, H.H. and Meuliet, E.J. Identification and development of mPGES-1 inhibitors: Where we are at? *Future Med. Chem.* **3(15)**, 1909-1934 (2011).
5. Korotkova, M. and Jakobsson, P.-J. Microsomal prostaglandin E synthase-1 in rheumatic diseases. *Frontiers in Pharmacology* **1(146)**, 1-8 (2011).

### 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, 12/02/2022

## 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