

# PRODUCT INFORMATION

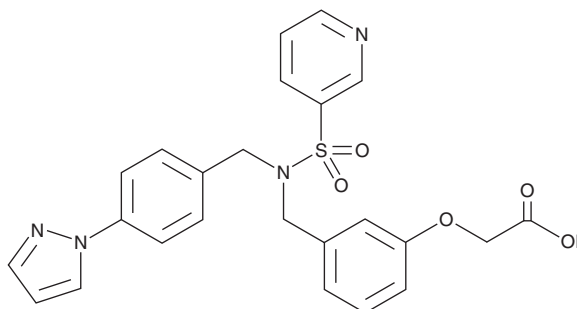


**CP 544,326**

Item No. 22945

**CAS Registry No.:** 752187-80-7  
**Formal Name:** 2-[3-[[[4-(1H-pyrazol-1-yl)phenyl]methyl](3-pyridinylsulfonyl)amino]methyl]phenoxy]-acetic acid

**MF:** C<sub>24</sub>H<sub>22</sub>N<sub>4</sub>O<sub>5</sub>S  
**FW:** 478.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 231, 262 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

CP 544,326 is supplied as a crystalline solid. A stock solution may be made by dissolving the CP 544,326 in the solvent of choice. CP 544,326 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of CP 544,326 in these solvents is approximately 30 and 20 mg/ml, respectively. CP 544,326 is also slightly soluble in ethanol.

## Description

CP 544,326 is a potent agonist of E prostanoid (EP) receptor 2 (IC<sub>50</sub>s = 10 and 15 nM for human and rat EP<sub>2</sub>, respectively).<sup>1</sup> It is selective for EP<sub>2</sub> over other EP receptors (IC<sub>50</sub>s = >3,200 nM for EP<sub>1</sub>, EP<sub>3</sub>, and EP<sub>4</sub>) and a panel of 37 G protein-coupled receptors (≤30% inhibition at a concentration of 10 μM). CP 544,326 increases cAMP levels in HEK293 cells expressing human EP<sub>2</sub> (EC<sub>50</sub> = 2.8 nM). CP 544,326 has poor corneal permeability in an *ex vivo* rabbit corneal model. However, PF-04217329, a cell-permeable prodrug form of CP 544,326, reduces intraocular pressure in rabbit, canine, and marmoset models of glaucoma.

## Reference

1. Prasanna, G., Carreiro, S., Anderson, S., *et al.* Effect of PF-04217329 a prodrug of a selective prostaglandin EP<sub>2</sub> agonist on intraocular pressure in preclinical models of glaucoma. *Exp. Eye Res.* **93(3)**, 256-264 (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, 12/22/2022

## 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