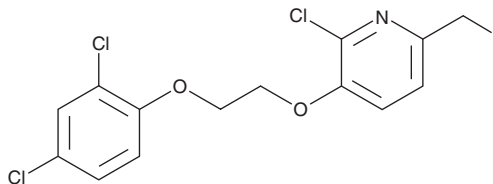


# PRODUCT INFORMATION



**CYM 50260**  
Item No. 14560

**CAS Registry No.:** 1355026-60-6  
**Formal Name:** 2-chloro-3-[2-(2,4-dichlorophenoxy)ethoxy]-6-(fluoromethyl)-pyridine  
**MF:** C<sub>14</sub>H<sub>11</sub>Cl<sub>3</sub>FNO<sub>2</sub>  
**FW:** 350.6  
**Purity:** ≥98%  
**Supplied as:** A 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

CYM 50260 is supplied as a solid. A stock solution may be made by dissolving the CYM 50260 in the solvent of choice, which should be purged with an inert gas. CYM 50260 is soluble (≥10 mg/ml) in DMSO.

## Description

CYM 50260 is an agonist of sphingosine-1-phosphate receptor 4 (S1P<sub>4</sub>).<sup>1</sup> It induces gene transcription in a reporter assay using U2OS cells expressing human S1P<sub>4</sub> (EC<sub>50</sub> = 45 nM). CYM 50260 (25 μM) inhibits collagen-induced increases in PDGF-AB secretion from, and phosphorylated heat shock protein 27 (Hsp27) levels in, isolated human platelets.<sup>2</sup> It also inhibits collagen-induced increases in platelet aggregate size.

## References

- Guerrero, M., Urbano, M., Zhao, J., *et al.* Discovery, design and synthesis of novel potent and selective sphingosine-1-phosphate 4 receptor (S1P<sub>4</sub>-R) agonists. *Bioorg. Med. Chem. Lett.* **22(1)**, 537-542 (2012).
- Onuma, T., Tanabe, K., Kito, Y., *et al.* Sphingosine 1-phosphate (S1P) suppresses the collagen-induced activation of human platelets via S1P4 receptor. *Thromb. Res.* **156**, 91-100 (2017).

### 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, 08/28/2025

## 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