

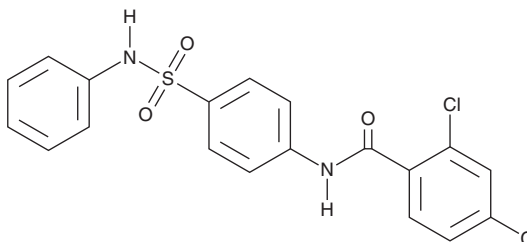
# PRODUCT INFORMATION



**CID-5157334**

Item No. 41677

**CAS Registry No.:** 651295-98-6  
**Formal Name:** 2,4-dichloro-N-[4-[(phenylamino) sulfonyl]phenyl]-benzamide  
**Synonym:** CID-5128535  
**MF:** C<sub>19</sub>H<sub>14</sub>Cl<sub>2</sub>N<sub>2</sub>O<sub>3</sub>S  
**FW:** 421.3  
**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

CID-5157334 is supplied as a solid. A stock solution may be made by dissolving the CID-5157334 in the solvent of choice, which should be purged with an inert gas. CID-5157334 is slightly soluble (0.1-1 mg/ml) in acetonitrile and sparingly soluble (1-10 mg/ml) in DMSO.

## Description

CID-5157334 is a surrogate agonist of G protein-coupled receptor 27 (GPR27).<sup>1</sup> It induces recruitment of  $\beta$ -arrestin-2 in complementation assays using wild-type GPR27 or a GPR27-vasopressin 2 (GPR27V<sub>2</sub>) chimeric receptor (EC<sub>50</sub>s = 7.943 and 0.417 nM, respectively) but does not induce recruitment of  $\beta$ -arrestin-1 in the same assay or activation of the G<sub>q</sub>, G<sub>s</sub>, or G<sub>i</sub> signaling pathways in aequorin-GFP assays. CID-5157334 is selective for GPR27 over GPR85V<sub>2</sub> and GPR173V<sub>2</sub> chimeric receptors at 15  $\mu$ M.

## Reference

1. Dupuis, N., Laschet, C., Franssen, D., *et al.* Activation of the orphan G protein-coupled receptor GPR27 by surrogate ligands promotes  $\beta$ -arrestin 2 recruitment. *Mol. Pharmacol.* **91**(6), 595-608 (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, 08/098/2024

## 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