

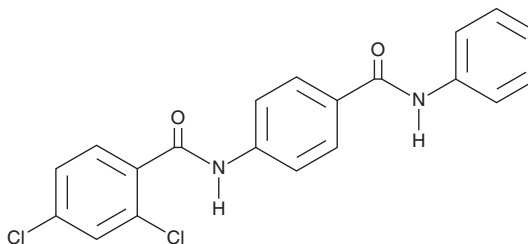
# PRODUCT INFORMATION



**CID-1375606**

Item No. 30422

**CAS Registry No.:** 313493-80-0  
**Formal Name:** 2,4-dichloro-N-[4-[(phenylamino) carbonyl]phenyl]-benzamide  
**MF:** C<sub>20</sub>H<sub>14</sub>Cl<sub>2</sub>N<sub>2</sub>O<sub>2</sub>  
**FW:** 385.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 290 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

## Laboratory Procedures

CID-1375606 is supplied as a crystalline solid. A stock solution may be made by dissolving the CID-1375606 in the solvent of choice, which should be purged with an inert gas. CID-1375606 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CID-1375606 in these solvents is approximately 5 and 10 mg/ml, respectively.

## Description

CID-1375606 is a GPR27 agonist with EC<sub>50</sub> values of 0.46 μM and 7.9 μM in chimeric GPR27V<sub>2</sub> and wild-type GPR27 luciferase β-arrestin-2 complementation assays, respectively.<sup>1</sup> It is selective for GPR27 over the closely related receptors GPR85 and GPR173 for which it does not induce β-arrestin-2 recruitment in the same assay. GPR27 function has been linked to insulin promoter activity and excretion.<sup>2</sup>

## References

1. Dupuis, N., Laschet, C., Franssen, D., *et al.* Activation of the orphan G protein-coupled receptor GPR27 by surrogate ligands promotes β-arrestin 2 recruitment. *Mol. Pharmacol.* **91(6)**, 595-608 (2017).
2. Ku, G.M., Pappalardo, Z., Luo, C.C., *et al.* An siRNA screen in pancreatic beta cells reveals a role for *Gpr27* in insulin production. *PLoS Genet.* **8(1)**, e1002449 (2012).

### 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, 05/26/2020

## 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