

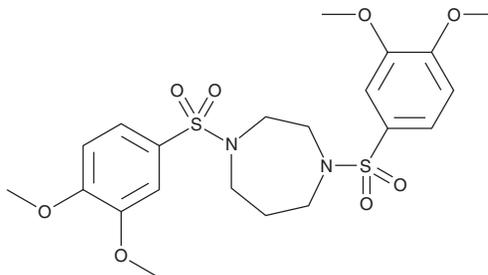
# PRODUCT INFORMATION



**SB-756050**

Item No. 27998

**CAS Registry No.:** 447410-57-3  
**Formal Name:** 1,4-bis[(3,4-dimethoxyphenyl)sulfonyl]hexahydro-1H-1,4-diazepine  
**MF:** C<sub>21</sub>H<sub>28</sub>N<sub>2</sub>O<sub>8</sub>S<sub>2</sub>  
**FW:** 500.6  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 249, 283 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

SB-756050 is supplied as a crystalline solid. A stock solution may be made by dissolving the SB-756050 in the solvent of choice, which should be purged with an inert gas. SB-756050 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of SB-756050 in these solvents is approximately 10 mg/ml.

SB-756050 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SB-756050 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. SB-756050 has a solubility of approximately 0.25 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

SB-756050 is an agonist of the bile acid G protein-coupled receptor TGR5 (EC<sub>50</sub> = 1.3 μM for the human receptor).<sup>1</sup> It is selective for TGR5 over the farnesoid X receptor (FXR) and a panel of other receptors, channels, and transporters. SB-756050 (10-100 mg/kg) decreases fasting glucose levels and increases glucose disposal rate and insulin secretion in diabetic Goto Kakizaki rats.

## Reference

1. Hodge, R.J., Lin, J., Vasist Johnson, L.S., *et al.* Safety, pharmacokinetics, and pharmacodynamic effects of a selective TGR5 agonist, SB-756050, in type 2 diabetes. *Clin. Pharmacol. Drug Dev.* **2**(3), 213-222 (2013).

### 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, 11/25/2019

## 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