

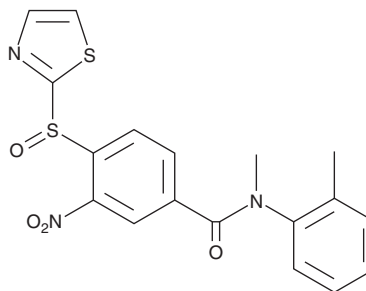
PRODUCT INFORMATION



SB-268262

Item No. 43547

CAS Registry No.: 217438-17-0
Formal Name: N-methyl-N-(2-methylphenyl)-3-nitro-4-(2-thiazolylsulfinyl)-benzamide
MF: C₁₈H₁₅N₃O₄S₂
FW: 401.5
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

SB-268262 is supplied as a solid. A stock solution may be made by dissolving the SB-268262 in the solvent of choice, which should be purged with an inert gas. SB-268262 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

SB-268262 is a nonpeptide calcitonin gene-related peptide (CGRP) receptor antagonist (IC₅₀ = 0.24 μM in a radioligand binding assay).¹ It inhibits CGRP-induced cAMP accumulation in SK-N-MC membrane preparations with an IC₅₀ value of 0.83 μM.

Reference

1. Aiyar, N., Daines, R.A., Disa, J., *et al.* Pharmacology of SB-273779, a nonpeptide calcitonin gene-related peptide 1 receptor antagonist. *J. Pharmacol. Exp. Ther.* **296**(3), 768-775 (2001).

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, 05/27/2025

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