

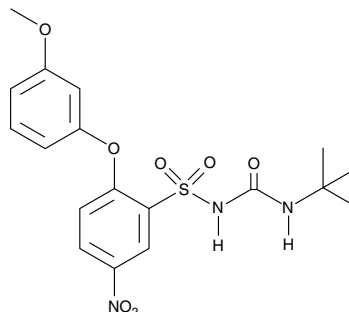
Product Information



CAY10535

Item No. 10010396

CAS Registry No.: 945716-28-9
Formal Name: N-(*tert*-butylcarbamoyl)-2-(3-methoxyphenoxy)-5-nitrobenzenesulfonamide
MF: C₁₈H₂₁N₃O₇S
FW: 423.4
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 294 nm



Laboratory Procedures

For long term storage, we suggest that CAY10535 be stored as supplied at -20°C. It should be stable for at least two years.

CAY10535 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10535 in an organic solvent purged with an inert gas. CAY10535 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CAY10535 in these solvents is approximately 10 mg/ml.

CAY10535 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10535 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CAY10535 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

TP α and TP β are two isoforms of the human TP receptor, the G protein-coupled receptor (GPCR) that mediates the actions of thromboxane A₂ (TXA₂). Although their distinct physiological functions have not been fully elucidated, TP β is believed to be responsible for vascular endothelial growth factor-induced endothelial cell differentiation and migration whereas TP α appears to be the predominant isoform expressed in platelets. CAY10535 is a TP receptor antagonist that shows ~20-fold selectivity for TP β (IC₅₀ = 99 nM) relative to TP α (IC₅₀ = 1,970 nM) in the inhibition of U-46619-mediated Ca²⁺ mobilization. This compound exhibits relatively poor activity on platelets (IC₅₀ = 985 nM) when inhibiting U-46619-induced platelet aggregation.¹

Reference

1. Hanson, J., Dogné, J.-M., Ghiotto, J., *et al.* Design, synthesis, and SAR study of a series of N-alkyl-N'-[2-(aryloxy)-5-nitrobenzenesulfonyl]ureas and -cyanoguanidine as selective antagonists of the TP α and TP β isoforms of the human thromboxane A₂ receptor. *J. Med. Chem.* **50**(16), 3928-3936 (2007).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10010396

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 10/03/2013

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com