

Product Information

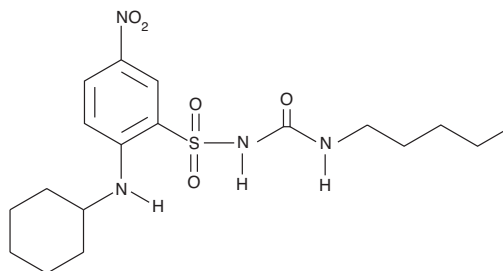


BM 567

Catalog No. 10155

Sold and made under non-exclusive license from Université de Liège.

CAS Registry No.: 284464-77-3
Formal Name: 2-(cyclohexylamino)-5-nitro-N-[(pentylamino)carbonyl]benzenesulfonamide
MF: C₁₈H₂₈N₄O₅S
FW: 412.5
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that BM 567 be stored as supplied at -20°C. It will be stable for at least two years.

BM 567 is supplied as a crystalline solid. A stock solution may be made by dissolving the BM 567 in an organic solvent purged with an inert gas. BM 567 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of BM 567 in these solvents is approximately 10 mg/ml in ethanol and 20 mg/ml in DMSO and DMF.

BM 567 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BM 567 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. BM 567 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Thromboxane A₂ (TXA₂) is a potent thrombogenic and vasoconstrictor eicosanoid, produced in large quantities by activated platelets. TXA₂ has been implicated as a causal factor in the onset of stroke and myocardial infarction. BM 567 is a dual acting antithrombogenic agent, acting as an inhibitor of TXA₂ synthase and as an antagonist of the TP receptor, the G protein-coupled receptor mediating TXA₂ activity in platelets and vascular smooth muscle. BM 567 antagonizes the vascular smooth muscle TP receptor with an IC₅₀ value of 1.1 nM. It inhibits platelet TX synthase with an IC₅₀ value of 12 nM.¹

Reference

1. Michaux, C., Rolin, S., Dogné, J.-M., *et al.* Structure determination and comparison of BM567, a sulfonylurea, with Terbogrel, two compounds with dual action, thromboxane receptor antagonism and thromboxane synthase inhibition. *Bioorg. Medicinal Chem. Letters* **11**, 1019-1022 (2001).

Related Products

BAY-u3405 - Cat. No. 10156 • U-44069 - Cat. No. 16440 • U-46619 - Cat. No. 16450 • U-51605 - Cat. No. 16465 • Pinane Thromboxane A₂ - Cat. No. 19020 • SQ 29,548 - Cat. No. 19025 • 1-Benzylimidazole - Cat. No. 70510 • Ozagrel - Cat. No. 70515 • Furegrelate (sodium salt) - Cat. No. 70540 • TP Receptor (murine) Polyclonal Antibody - Cat. No. 101882 • Thromboxane Synthase Polyclonal Antibody - Cat. No. 160715 • TP Receptor (human) Polyclonal Antibody - Cat. No. 10004452

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

WARNING: THIS PRODUCT IS NOT INTENDED OR APPROVED FOR HUMAN OR VETERINARY USE. USE OF THIS PRODUCT FOR HUMAN OR ANIMAL TESTING IS EXTREMELY HAZARDOUS AND MAY RESULT IN DISEASE, SEVERE INJURY, OR DEATH.

MATERIAL 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 Material Safety Data Sheet, which has been sent under separate cover to the MSDS supervisor at 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, 03/29/2007