

PRODUCT INFORMATION

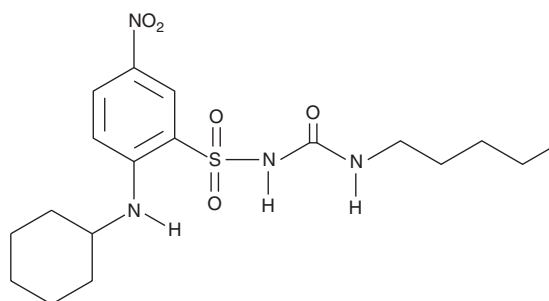


BM 567

Item No. 10155

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

CAS Registry No.: 284464-77-3
Formal Name: 2-(cyclohexylamino)-5-nitro-N-[(pentylamino)carbonyl]benzenesulfonamide
MF: C₁₈H₂₈N₄O₅S
FW: 412.5
Purity: ≥98%
UV/Vis.: λ_{max}: 212, 365 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

BM 567 is supplied as a crystalline solid. A stock solution may be made by dissolving the BM 567 in the solvent of choice, which should be 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 ethanol is approximately 10 mg/ml and approximately 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.

Description

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. Med. Chem. Lett.* **11(8)**, 1019-1022 (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, 11/07/2023

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