

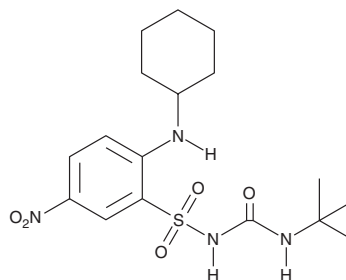
PRODUCT INFORMATION



BM-531

Item No. 45368

CAS Registry No.: 284464-46-6
Formal Name: 2-(cyclohexylamino)-N-[[[(1,1-dimethylethyl)amino]carbonyl]-5-nitro-benzenesulfonamide
MF: C₁₇H₂₆N₄O₅S
FW: 398.5
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

BM-531 is supplied as a solid. A stock solution may be made by dissolving the BM-531 in the solvent of choice, which should be purged with an inert gas. BM-531 is slightly soluble (0.1-1 mg/ml) in ethanol and sparingly soluble (1-10 mg/ml) in DMSO.

Description

BM-531 is an antagonist of the thromboxane A₂ (TXA₂) receptor (TP receptor; IC₅₀ = 0.0078 μM) and an inhibitor of thromboxane synthase.^{1,2} It inhibits the production of TXB₂ by arachidonic acid-activated human platelets when used at concentrations of 1 or 10 μM.² BM-531 prevents platelet aggregation induced by arachidonic acid (Item No. 90010; ED₁₀₀ = 0.125 μM) and U-46619 (Item No. 16450; ED₅₀ = 0.482 μM) in human citrated platelet-rich plasma.¹ It also prevents collagen-induced platelet aggregation by 42.9% in human citrated platelet-rich plasma when used at a concentration of 10 μM. BM-531 reduces U-46619-induced fundus contractions in rats (ED₅₀ = 1.5 μM) but not spontaneous contractions or those induced by prostaglandin E₂ (PGE₂; Item No. 14010), PGF_{2α} (Item Nos. 16010 | 16020), or PGI₂ (Item No. 18220).²

References

1. Dogné, J.M., de Leval, X., Kolh, P., *et al.* Pharmacological evaluation of the novel thromboxane modulator BM-567 (I/II). Effects of BM-567 on platelet function. *Prostaglandins Leukot. Essent. Fatty Acids* **68**(1), 49-54 (2003).
2. Dogné, J.M., Rolin, S., de Leval, X., *et al.* Pharmacology of the thromboxane receptor antagonist and thromboxane synthase inhibitor BM-531. *Cardiovasc. Drug Rev.* **19**(2), 87-96 (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, 04/03/2026

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