

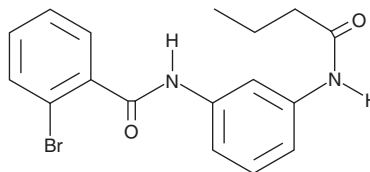
PRODUCT INFORMATION



ML-161

Item No. 15179

CAS Registry No.: 423735-93-7
Formal Name: 2-bromo-N-[3-[(1-oxobutyl)amino]phenyl]-benzamide
Synonyms: AG-670, CID-1048267
MF: C₁₇H₁₇BrN₂O₂
FW: 361.2
Purity: ≥95%
UV/Vis.: λ_{max}: 240 nm
Supplied as: A crystalline 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

ML-161 is supplied as a crystalline solid. A stock solution may be made by dissolving the ML-161 in the solvent of choice. ML-161 is soluble in organic solvents such as ethanol and DMSO, which should be purged with an inert gas. The solubility of ML-161 in these solvents is approximately 25 mg/ml.

Description

ML-161 is an allosteric, reversible inhibitor of proteinase-activated receptor 1 (PAR1) on platelets, preventing surface expression of P-selectin induced by the peptide SFLLRN with an IC₅₀ value of 0.26 μM.¹ It blocks platelet activation induced by thrombin as well as by SFLLRN but not by PMA (Item No. 10008014), U-46619 (Item No. 16450), or collagen.¹

Reference

1. Dockendorff, C., Aisiku, O., VerPlank, L., *et al.* Discovery of 1,3-diaminobenzenes as selective inhibitors of platelet activation at the PAR1 receptor. *ACS Med. Chem. Lett.* **3**(3), 232-237 (2012).

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, 10/04/2022

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