

PRODUCT INFORMATION

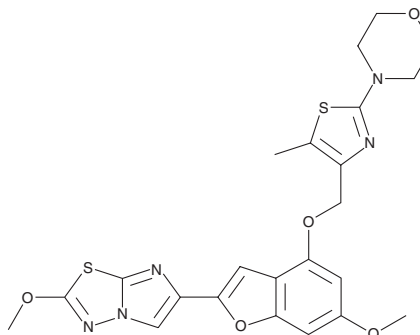


BMS 986120

Item No. 23497

CAS Registry No.: 1478712-37-6
Formal Name: 2-methoxy-6-[6-methoxy-4-[[5-methyl-2-(4-morpholinyl)-4-thiazolyl]methoxy]-2-benzofuranyl]-imidazo[2,1-b]-1,3,4-thiadiazole

MF: C₂₃H₂₃N₅O₅S₂
FW: 513.6
Purity: ≥98%
UV/Vis.: λ_{max}: 305 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

BMS 986120 is supplied as a crystalline solid. A stock solution may be made by dissolving the BMS 986120 in the solvent of choice, which should be purged with an inert gas. BMS 986120 is soluble in the organic solvent DMSO.

Description

BMS 986120 is an orally bioavailable, selective, and reversible antagonist of proteinase-activated receptor 4 (PAR4; IC₅₀ = 0.56 nM to inhibit calcium mobilization induced by PAR4 agonist peptide (PAR4-AP) in HEK293 cells).¹ It is selective for PAR4 over PAR1, PAR2, and a panel of purified proteases, including thrombin. It inhibits platelet aggregation *in vitro* in human platelet-rich plasma (IC₅₀ = 7.3 nM). BMS 986120 (0.2-1 mg/kg) decreases platelet aggregation induced by PAR4-AP *ex vivo* in a dose-dependent manner, but does not increase clotting times. *In vivo*, BMS 986120 (1 mg/kg) prevents vascular occlusion and reduces thrombus formation by 82% in cynomolgus monkeys.

Reference

1. Wong, P.C., Seiffert, D., Bird, J.E., *et al.* Blockade of protease-activated receptor-4 (PAR4) provides robust antithrombotic activity with low bleeding. *Sci. Transl. Med.* **9**(371), eaaf5294 (2017).

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, 10/26/2022

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