

PRODUCT INFORMATION

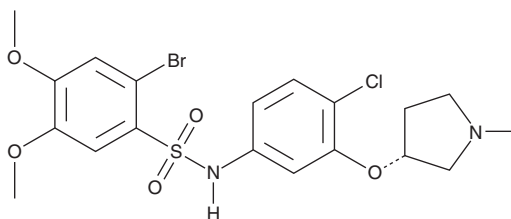


SB-657510

Item No. 42321

CAS Registry No.: 474960-44-6
Formal Name: 2-bromo-N-[[4-chloro-3-[[[(3R)-1-methyl-3-pyrrolidinyl]oxy]phenyl]-4,5-dimethoxybenzenesulfonamide

MF: C₁₉H₂₂BrClN₂O₅S
FW: 505.8
Purity: ≥95%
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

SB-657510 is supplied as a solid. A stock solution may be made by dissolving the SB-657510 in the solvent of choice, which should be purged with an inert gas. SB-657510 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

SB-657510 is an antagonist of the urotensin II receptor (UTR; K_is = 61, 17, 30, 65, and 56 nM for the human, monkey, cat, rat, and mouse receptors, respectively).¹ It reduces urotensin II-induced increases in the expression of *IL1B*, *IL6*, and *VCAM1* in EA.hy926 human endothelial cells and urotensin II-induced adhesion of U937 human monocytes to EA.hy926 cells when used at a concentration of 1 μM.²

References

1. Behm, D.J., McAtee, J.J., Dodson, J.W., *et al.* Palosuran inhibits binding to primate UT receptors in cell membranes but demonstrates differential activity in intact cells and vascular tissues. *Brit. J. Pharmacol.* **155**(3), 374-386 (2008).
2. Park, S.L., Lee, B.K., Kim, Y.-A., *et al.* Inhibitory effect of an urotensin II receptor antagonist on proinflammatory activation induced by urotensin II in human vascular endothelial cells. *Biomol. Ther. (Seoul)* **21**(4), 277-283 (2013).

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, 11/11/2024

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