

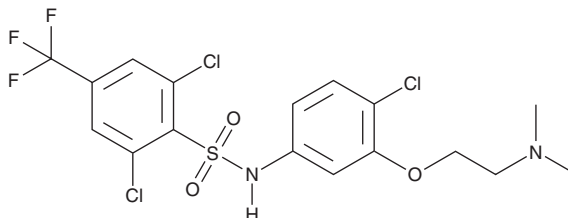
PRODUCT INFORMATION



SB-611812

Item No. 33247

CAS Registry No.: 345892-71-9
Formal Name: 2,6-dichloro-N-[4-chloro-3-[2-(dimethylamino)ethoxy]phenyl]-4-(trifluoromethyl)benzenesulfonamide
MF: C₁₇H₁₆Cl₃F₃N₂O₃S
FW: 491.7
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

SB-611812 is supplied as a solid. A stock solution may be made by dissolving the SB-611812 in the solvent of choice, which should be purged with an inert gas. SB-611812 is soluble in the organic solvent DMSO at a concentration of approximately 10 mM.

Description

SB-611812 is an antagonist of the urotensin II receptor (UTR).¹ It binds to recombinant rat UTR ($K_i = 121$ nM) and inhibits contraction of isolated rat aorta induced by urotensin II ($pA_2 = 6.59$) but not potassium chloride, endothelin-1, angiotensin II, or phenylephrine. SB-611812 (30 mg/kg) decreases myocardial interstitial fibrosis and the ratio of collagen I to collagen III in a rat model of ischemia induced by ligation of the left anterior descending (LAD) artery.²

References

1. Rakowski, E., Hassan, G.S., Dhanak, D., *et al.* A role for urotensin II in restenosis following balloon angioplasty: Use of a selective UT receptor blocker. *J. Mol. Cell. Cardiol.* **39(5)**, 785-791 (2005).
2. Bousette, N., Pottinger, J., Ramli, W., *et al.* Urotensin-II receptor blockade with SB-611812 attenuates cardiac remodeling in experimental ischemic heart disease. *Peptides* **27(11)**, 2919-2926 (2006).

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, 11/14/2022

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