

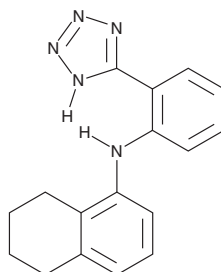
PRODUCT INFORMATION



BL 1249

Item No. 37426

CAS Registry No.: 18200-13-0
Formal Name: 5,6,7,8-tetrahydro-N-[2-(2H-tetrazol-5-yl)phenyl]-1-naphthalenamine
MF: C₁₇H₁₇N₅
FW: 291.4
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

BL 1249 is supplied as a solid. A stock solution may be made by dissolving the BL 1249 in the solvent of choice, which should be purged with an inert gas. BL 1249 is slightly soluble (0.1-1 mg/ml) in acetonitrile and methanol, and sparingly soluble (1-10 mg/ml) in DMSO.

Description

BL 1249 is an activator of mechanosensitive two-pore domain potassium channels (K_{2p}).¹ It selectively induces currents in *Xenopus* oocytes expressing K_{2p}2.1/TREK-1 or K_{2p}10.1/TREK-2 over *Xenopus* oocytes expressing K_{2p}4.1/TRAAK, K_{2p}3.1/TASK1, K_{2p}3.2/TASK2, K_{2p}3.3/TASK3, K_{2p}18.1/TRESK, K_{2p}13.1/THIK1, or K_{2p}1.1/TWIK-1 at 10 μM.

Reference

1. Pope, L., Arrigoni, C., Lou, H., *et al.* Protein and chemical determinants of BL-1249 action and selectivity for K_{2p} channels. *ACS Chem. Neurosci.* **9**(12), 3153-3165 (2018).

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, 06/04/2024

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