

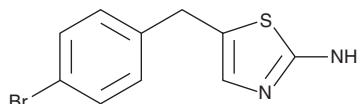
PRODUCT INFORMATION



NS 19504

Item No. 43669

CAS Registry No.: 327062-46-4
Formal Name: 5-[(4-bromophenyl)methyl]-2-thiazolamine
MF: C₁₀H₉BrN₂S
FW: 269.2
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

NS 19504 is supplied as a solid. A stock solution may be made by dissolving the NS 19504 in the solvent of choice, which should be purged with an inert gas. NS 19504 is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

NS 19504 is an activator of large-conductance calcium-activated potassium channels (K_{Ca}1.1/BK).¹ It induces currents in whole-cell patch-clamp assays using HEK293 cells expressing human K_{Ca}1.1/BK when used at a concentration of 10 μM. NS 19504 decreases the force of spontaneous phasic contractions in isolated and perfused guinea pig bladder smooth muscle strips (EC₅₀ = 640 nM). It inhibits methoxamine-induced contractions in isolated rat saphenous arteries when used at a concentration of 6 μM.²

References

1. Nausch, B., Rode, F., Jørgensen, S., *et al.* NS19504: A novel BK channel activator with relaxing effect on bladder smooth muscle spontaneous phasic contractions. *J. Pharmacol. Exp. Ther.* **350(3)**, 520-530 (2014).
2. Ma, D., Gaynullina, D., Schmidt, N., *et al.* The functional availability of arterial Kv7 channels is suppressed considerably by large-conductance calcium-activated potassium channels in 2- to 3-month old but not in 10- to 15-day old rats. *Front. Physiol.* **11**, 597395 (2020).

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, 08/04/2025

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