

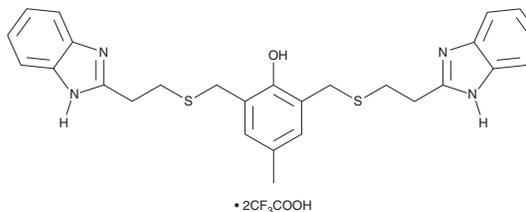
PRODUCT INFORMATION



NSC 620023 (trifluoroacetate salt)

Item No. 401037

Formal Name: 2,6-bis[[[2-(1H-benzimidazol-2-yl)ethyl]thio]methyl]-4-methylphenol, trifluoroacetate salt
MF: C₂₇H₂₈N₄OS₂ • 2CF₃COOH
FW: 716.7
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

NSC 620023 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the NSC 620023 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. NSC 620023 (trifluoroacetate salt) is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

NSC 620023 is an inhibitor of butyrylcholinesterase (BChE; IC₅₀ = 40 nM).¹ It is selective for BChE over acetylcholinesterase (AChE; IC₅₀ = 2,604 nM).

Reference

1. Williams, A., Zhou, S., and Zhan, C.G. Discovery of potent and selective butyrylcholinesterase inhibitors through the use of pharmacophore-based screening. *Bioorg. Med. Chem. Lett.* **29(24)**, 126754 (2019).

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, 07/31/2025

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