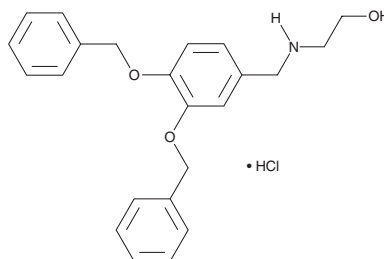


PRODUCT INFORMATION



XIE62-1004
Item No. 40446

CAS Registry No.: 2421146-32-7
Formal Name: 2-[[[3,4-bis(phenylmethoxy)phenyl]methyl]amino]-ethanol, monohydrochloride
MF: C₂₃H₂₅NO₃ • HCl
FW: 399.9
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

XIE62-1004 is supplied as a solid. A stock solution may be made by dissolving the XIE62-1004 in the solvent of choice, which should be purged with an inert gas. XIE62-1004 is soluble in DMSO. XIE62-1004 is slightly soluble in acetonitrile.

XIE62-1004 is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

XIE62-1004 is an inducer of autophagy.¹ It increases the number of p62 aggregates, as well as the number of autophagosomes and autolysosomes, in wild-type HeLa cervical cancer cells, but not in HeLa cells deficient in the gene encoding p62, *SQSTM1*, when used at a concentration of 10 μM. XIE62-1004 induces autophagic flux and the conversion of LC3 autophagy marker I (LC3-I) to LC3-II in HeLa cells. It increases the degradation of mutant huntingtin (mHTT^{Q103}) protein expressed in HeLa cells.

Reference

1. Cha-Molstad, H., Yu, J.E., Feng, Z., *et al.* p62/SQSTM1/Sequestosome-1 is an N-recogin of the N-end rule pathway which modulates autophagosome biogenesis. *Nat. Commun.* **8(1)**, 102 (2017).

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, 03/05/2024

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