

PRODUCT INFORMATION

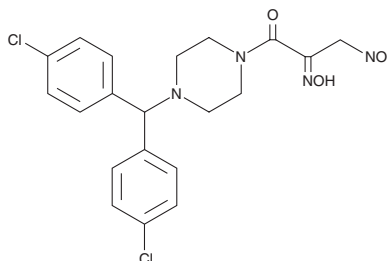


JKE-1674

Item No. 30784

CAS Registry No.: 2421119-60-8
Formal Name: (E)-1-(4-(bis(4-chlorophenyl)methyl)piperazin-1-yl)-2-(hydroxyimino)-3-nitropropan-1-one

MF: C₂₀H₂₀Cl₂N₄O₄
FW: 451.3
Purity: ≥98% (mixture of isomers)
UV/Vis.: λ_{max}: 232 nm
Supplied as: A crystalline 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

JKE-1674 is supplied as a crystalline solid. A stock solution may be made by dissolving the JKE-1674 in the solvent of choice, which should be purged with an inert gas. JKE-1674 is soluble in DMSO.

Description

JKE-1674 is an inhibitor of glutathione peroxidase 4 (GPX4) and an active metabolite of the GPX4 inhibitor ML-210 (Item No. 23282).¹ JKE-1674 reduces viability of LOX-IMVI cancer cells (EC₅₀ = 0.03 μM) and in a panel of additional cancer cell lines, an effect that can be blocked by the ferroptosis inhibitor ferrostatin-1 (Item No. 17729).

Reference

1. Eaton, J.K., Furst, L., Ruberto, R.A., *et al.* Selective covalent targeting of GPX4 using masked nitrile-oxide electrophiles. *Nat. Chem. Biol.* **16**(5), 497-506 (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, 03/21/2024

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