

# PRODUCT INFORMATION



## JKE-1716

Item No. 30786

CAS Registry No.: 2421118-05-8

Formal Name: 3-[4-[bis(4-chlorophenyl)methyl]-1-piperazinyl]-1-nitro-1,3-propanedione, 1-oxime

MF: C<sub>20</sub>H<sub>20</sub>Cl<sub>2</sub>N<sub>4</sub>O<sub>4</sub>

FW: 451.3

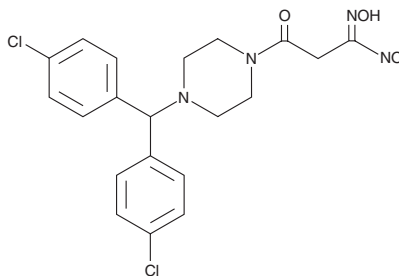
Purity: ≥98%

UV/Vis.: λ<sub>max</sub>: 233 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

JKE-1716 is supplied as a crystalline solid. A stock solution may be made by dissolving the JKE-1716 in the solvent of choice, which should be purged with an inert gas. JKE-1716 is soluble in DMSO.

### Description

JKE-1716 is an inhibitor of glutathione peroxidase 4 (GPX4) and a derivative of the GPX4 inhibitor ML-210 (Item No. 23282).<sup>1</sup> JKE-1716 reduces viability of LOX-IMVI cancer cells in a concentration-dependent manner 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, 11/06/2020

#### 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