

PRODUCT INFORMATION



PKUMDL-LC-101-D04

Item No. 28966

CAS Registry No.: 2143896-83-5
Formal Name: N-(2-aminoethyl)-4-[[[(cyclopentylamino)thioxomethyl]amino]-benzenesulfonamide, monohydrochloride

MF: C₁₄H₂₂N₄O₂S₂ • HCl

FW: 378.9

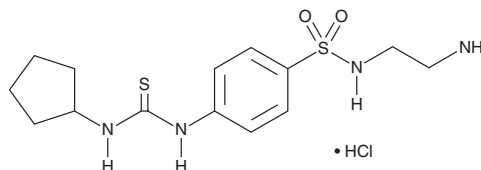
Purity: ≥98%

UV/Vis.: λ_{max}: 260, 291 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

PKUMDL-LC-101-D04 is supplied as a crystalline solid. A stock solution may be made by dissolving the PKUMDL-LC-101-D04 in the solvent of choice, which should be purged with an inert gas. PKUMDL-LC-101-D04 is soluble in the organic solvent DMSO.

Description

PKUMDL-LC-101-D04 is an allosteric activator of glutathione peroxidase 4 (GPX4).¹ It increases GPX4 activity to 150% of control levels when used at a concentration of 20 μM in a cell-free assay and at 61 μM in wild-type, but not *Gpx4*^{-/-}, mouse embryonic fibroblast (MEF) extracts. PKUMDL-LC-101-D04 (200 μM) also reduces cholesterol hydroperoxide-induced MEF death.

Reference

1. Li, C., Deng, X., Zhang, W., *et al.* Novel allosteric activators for ferroptosis regulator glutathione peroxidase 4. *J. Med. Chem.* **62**(1), 266-275 (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, 10/19/2022

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