

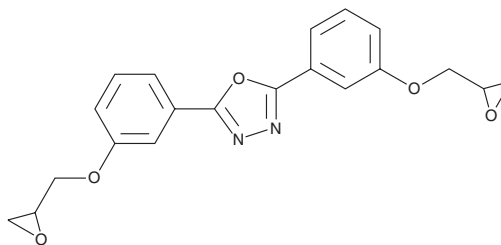
PRODUCT INFORMATION



GAPDH Inhibitor Compound F8

Item No. 42090

CAS Registry No.: 252212-58-1
Formal Name: 2,5-bis[3-(2-oxiranylmethoxy)phenyl]-1,3,4-oxadiazole
MF: C₂₀H₁₈N₂O₅
FW: 366.4
Purity: ≥95%
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

GAPDH inhibitor compound F8 is supplied as a solid. A stock solution may be made by dissolving the GAPDH inhibitor compound F8 in the solvent of choice, which should be purged with an inert gas. GAPDH inhibitor compound F8 is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

GAPDH inhibitor compound F8 is a covalent inhibitor of GAPDH (IC₅₀ = 39.31 μM).¹ It decreases viability of, and glucose metabolism in, HEK293 cells in a concentration-dependent manner.

Reference

1. Jung, Y., Noda, N., Takaya, J., *et al.* Discovery of non-cysteine-targeting covalent inhibitors by activity-based proteomic screening with a cysteine-reactive probe. *ACS Chem. Biol.* **17(2)**, 340-347 (2022).

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, 10/28/2024

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