

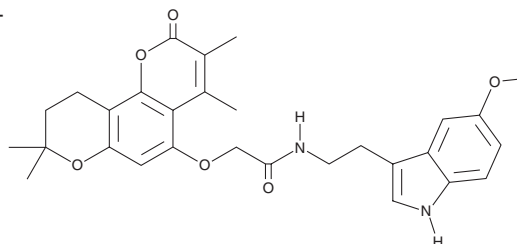
PRODUCT INFORMATION



NPD4928

Item No. 36598

CAS Registry No.: 956965-61-0
Formal Name: 2-[(9,10-dihydro-3,4,8,8-tetramethyl-2-oxo-2H,8H-benzo[1,2-b:3,4-b']dipyran-5-yl)oxy]-N-[2-(5-methoxy-1H-indol-3-yl)ethyl]-acetamide
MF: C₂₉H₃₂N₂O₆
FW: 504.6
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

NPD4928 is supplied as a solid. A stock solution may be made by dissolving the NPD4928 in the solvent of choice, which should be purged with an inert gas. NPD4928 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

NPD4928 is an inhibitor of ferroptosis suppressor protein 1 (FSP1) and an enhancer of ferroptosis.¹ It binds to FSP1 and inhibits its activity in a cell-free NADH consumption assay. NPD4928 (3 μM) enhances cytotoxicity induced by the ferroptosis inducer RSL3 (Item No. 19288) in PANC-1 prostate cancer cells in a synergistic manner. It also enhances cytotoxicity induced by sulfasalazine (Item No. 15025) and L-buthionine sulfoximine (BSO) in the same cells. NPD4928 in combination with RSL3 induces cytotoxicity in PC3, MKN74, DU145, MIA PaCa-2, and A549 cancer cells (IC₅₀s = 105, 95, 103, 140, and 155 nM, respectively) but not BxPC-3 or RPMI-8226 cancer cells (IC₅₀s = >30 μM for both).

Reference

1. Yoshioka, H., Kawamura, T., Muroi, M., *et al.* Identification of a small molecule that enhances ferroptosis via inhibition of ferroptosis suppressor protein 1 (FSP1). *ACS Chem Biol.* **17**(2), 483-491 (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, 11/05/2024

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