

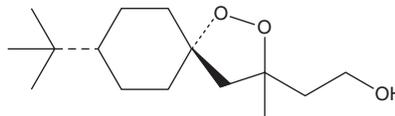
PRODUCT INFORMATION



FINO₂

Item No. 25096

CAS Registry No.: 869298-31-7
Formal Name: (5 α ,8 α)-8-(1,1-dimethylethyl)-3-methyl-1,2-dioxaspiro[4.5]decane-3-ethanol
MF: C₁₅H₂₈O₃
FW: 256.4
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

FINO₂ is supplied as a solid. A stock solution may be made by dissolving the FINO₂ in the solvent of choice. FINO₂ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of FINO₂ in these solvents is approximately 30 mg/ml.

FINO₂ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, FINO₂ should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. FINO₂ has a solubility of approximately 0.3 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

FINO₂ is a ferroptosis-inducing peroxide compound that indirectly inhibits glutathione peroxidase 4 (GPX4) and oxidizes iron.¹ It decreases GPX4 activity and protein levels *in vitro* but does not act as an active site, allosteric, or covalent inhibitor of GPX4 or alter GPX homeostasis. It also oxidizes iron *in vitro*, leading to degradation of the endoperoxide moiety, but does not affect the protein levels of iron regulatory proteins, such as IRP2, FTL1, or TFR. FINO₂ induces lipid peroxidation of a large subset of the lipidome in HT-1080 cells when used at a concentration of 10 μ M and induces ferroptosis in an arachidonic acid lipoxygenase-independent manner. It inhibits cell growth (mean GI₅₀ = 5.8 μ M) and induces lethality (mean LC₅₀ = 46 μ M) in the NCI-60 panel of cancer cell lines.² It is selective for oncogenically transformed BJ-ELR cells over noncancerous BJ-hTERT cells when used at concentrations of 15 and 20 μ M. FINO₂ (6 μ M) induces oxidative stress, including lipid peroxidation, in RS4;11 B-lymphoblastic leukemia cells. It induces iron-dependent cell death, an effect that can be blocked by pretreatment with the lipophilic antioxidants ferrostatin-1 (Item No. 17729) and liproxstatin-1 (Item No. 17730), and does not induce markers of apoptosis, necrosis, or autophagy in RS4;11 cells.

References

1. Gaschler, M.M., Andia, A.A., Liu, H., *et al.* FINO₂ initiates ferroptosis through GPX4 inactivation and iron oxidation. *Nat. Chem. Biol.* **14**(5), 507-515 (2018).
2. Abrams, R.P., Carroll, W.L., and Woerpel, K.A. Five-membered ring peroxide selectively initiates ferroptosis in cancer cells. *ACS Chem. Biol.* **11**(5), 1305-1312 (2016).

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, 09/28/2022

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