

# PRODUCT INFORMATION



## Ferroptocide

Item No. 39005

**CAS Registry No.:** 2505218-38-0  
**Formal Name:** (2S,3aS,4S,5R,7R,8S,8aS,9R)-octahydro-2,8a-dihydroxy-4,8,9-trimethyl-1-oxo-5-(2,3,5,8-tetrahydro-1,3-dioxo-2-phenyl-1H-[1,2,4]triazolo[1,2-a]pyridazin-6-yl)-1H-3a,8-propanoazulen-7-yl ester, 2-chloro-acetic acid

**MF:** C<sub>30</sub>H<sub>36</sub>ClN<sub>3</sub>O<sub>7</sub>

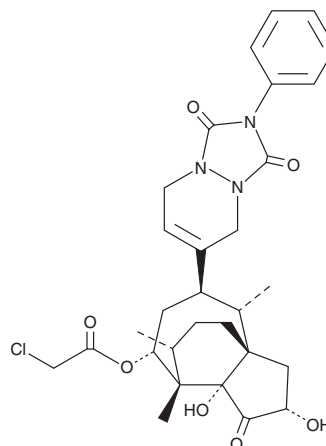
**FW:** 586.1

**Purity:** ≥90%

**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

Ferroptocide is supplied as a solid. A stock solution may be made by dissolving the ferroptocide in the solvent of choice, which should be purged with an inert gas. Ferroptocide is soluble in DMSO and methanol.

### Description

Ferroptocide is an inhibitor of thioredoxin (Trx) and a derivative of pleuromutilin (Item No. 19452).<sup>1</sup> It inhibits Trx activity in ES-2 ovarian cancer cells when used at a concentration of 20 μM and is cytotoxic to ES-2 cells (IC<sub>50</sub> = 1.6 μM). Ferroptocide (5 μM) induces ferroptosis in ES-2 cells, an effect that can be blocked by the iron chelator deferoxamine (DFO; Item No. 14595), antioxidant Trolox (Item No. 10011659), or ferroptosis inhibitor ferrostatin-1 (Item No. 17729). It decreases tumor volume in a 4T1 murine mammary carcinoma model using immunocompetent but not immunodeficient mice when administered at 50 mg/kg twice per week.

### Reference

1. Llabani, E., Hicklin, R.W., Lee, H.Y., *et al.* Diverse compounds from pleuromutilin lead to a thioredoxin inhibitor and inducer of ferroptosis. *Nat. Chem.* **11**(6), 521-532 (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, 10/19/2023

#### 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