

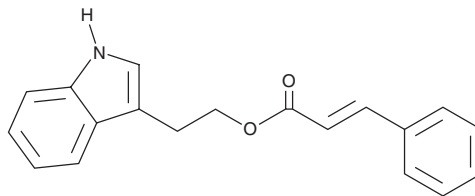
# PRODUCT INFORMATION



## Ferroptosis Inducer 24

Item No. 41201

CAS Registry No.: 28657-88-7  
Formal Name: 3-phenyl-2-propenoic acid, 2-(1H-indol-3-yl)ethyl ester  
MF:  $C_{19}H_{17}NO_2$   
FW: 291.3  
Purity:  $\geq 95\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

Ferroptosis inducer 24 is supplied as a solid. A stock solution may be made by dissolving the ferroptosis inducer 24 in the solvent of choice, which should be purged with an inert gas. Ferroptosis inducer 24 is soluble ( $\geq 10$  mg/ml) in DMSO and sparingly soluble (1-10 mg/ml) in ethanol.

### Description

Ferroptosis inducer 24 is an inducer of ferroptosis.<sup>1</sup> It inhibits glutathione peroxidase 4 (GPX4) by 30-35% when used at a concentration of 50  $\mu\text{M}$ . Ferroptosis inducer 24 (10  $\mu\text{M}$ ) increases the levels of heme oxygenase-1 (HO-1) in MDA-MB-231 triple-negative breast cancer cells, as well as increases the levels of bilirubin formed by HO-1 in MDA-MB-231 cell lysates. It reduces the viability of, as well as increases the levels of oxidized lipids, decreases the levels of IL-6 and glutathione (GST), and disrupts the mitochondrial membrane potential in, MDA-MB-231 cancer cells when used at a concentration of 50  $\mu\text{M}$ . Ferroptosis inducer 24 also increases the levels of reactive oxygen species (ROS) in a time-dependent manner.

### Reference

1. Consoli, V., Fallica, A.N., Virzi, N.F., *et al.* Synthesis and in vitro evaluation of CAPE derivatives as ferroptosis inducers in triple negative breast cancer. *ACS Med. Chem. Lett.* **15**(5), 706-713 (2024).

#### 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, 06/10/2024

#### 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