

# PRODUCT INFORMATION



## viFSP1

Item No. 39927

**CAS Registry No.:** 951945-67-8  
**Formal Name:** 5-amino-1-(2,4-dimethoxyphenyl)-1,2-dihydro-4-(4-methyl-2-thiazolyl)-3H-pyrrol-3-one

**Synonyms:** Versatile Inhibitor of Ferroptosis Suppressor Protein 1, Versatile Inhibitor of FSP1

**MF:** C<sub>16</sub>H<sub>17</sub>N<sub>3</sub>O<sub>3</sub>S

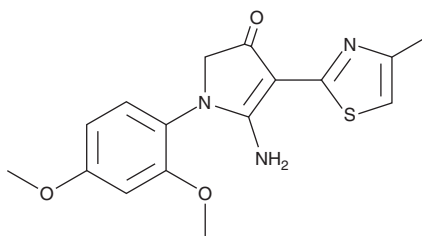
**FW:** 331.4

**Purity:** ≥95%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥3 years



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

### Laboratory Procedures

viFSP1 is supplied as a solid. A stock solution may be made by dissolving the viFSP1 in the solvent of choice, which should be purged with an inert gas. viFSP1 is soluble in acetonitrile and chloroform.

### Description

viFSP1 is a species-independent inhibitor of ferroptosis suppressor protein 1 (FSP1; IC<sub>50</sub>s = 34 and 83 nM for the human and mouse enzymes, respectively).<sup>1</sup> It induces ferroptosis in glutathione peroxidase 4 (Gpx4) knockout Pfa1 mouse embryonic fibroblasts expressing human or mouse FSP1, an effect that can be reversed by the ferroptosis inhibitor liproxstatin-1 (Item No. 17730).

### Reference

1. Nakamura, T., Mishima, E., Yamada, N., *et al.* Integrated chemical and genetic screens unveil FSP1 mechanisms of ferroptosis regulation. *Nat. Struct. Mol. Biol.* **30(11)**, 1806-1815 (2023).

#### 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, 02/21/2024

#### 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