

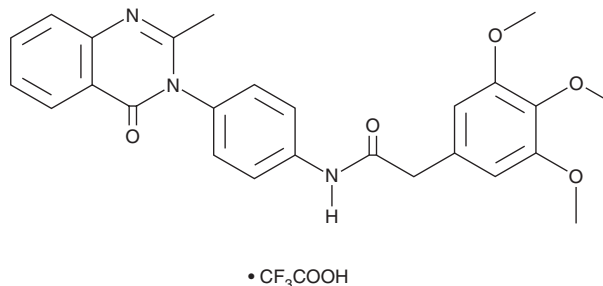
# PRODUCT INFORMATION



## icFSP1 (trifluoroacetate salt)

Item No. 39347

**Formal Name:** 3,4,5-trimethoxy-N-[4-(2-methyl-4-oxo-3(4H)-quinazoliny)phenyl]-benzeneacetamide, trifluoroacetate salt  
**MF:** C<sub>26</sub>H<sub>25</sub>N<sub>3</sub>O<sub>5</sub> • CF<sub>3</sub>COOH  
**FW:** 573.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 236 nm  
**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.

### Description

icFSP1 is an inhibitor of ferroptosis suppressor protein 1 (FSP1).<sup>1</sup> It induces condensation and phase separation of FSP1 in cells when used at a concentration of 2.5 μM but does not inhibit FSP1 enzymatic activity in a cell-free assay (IC<sub>50</sub> = > 30 μM). icFSP1 induces ferroptosis in HT-1080 fibrosarcoma cells in a concentration-dependent manner. *In vivo*, icFSP1 (50 mg/kg) decreases tumor volume and weight in a B16/F10 murine melanoma model using human FSP1-overexpressing, *Gpx4*<sup>-/-</sup>*Fsp1*<sup>-/-</sup> B16/F10 cells.

### Reference

1. Nakamura, T., Hipp, C., Mourão, A.S.D., et al. Phase separation of FSP1 promotes ferroptosis. *Nature* **619**, 371-377 (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, 09/20/2023

#### 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