

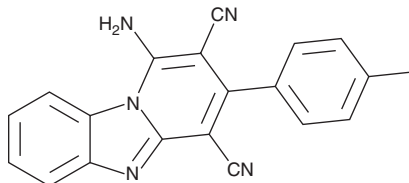
PRODUCT INFORMATION



iFSP1

Item No. 29483

CAS Registry No.: 150651-39-1
Formal Name: 1-amino-3-(4-methylphenyl)-pyrido[1,2-a]benzimidazole-2,4-dicarbonitrile
MF: C₂₀H₁₃N₅
FW: 323.4
Purity: ≥98%
UV/Vis.: λ_{max}: 248, 283 nm
Supplied as: A crystalline 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

iFSP1 is supplied as a crystalline solid. A stock solution may be made by dissolving the iFSP1 in the solvent of choice, which should be purged with an inert gas. iFSP1 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of iFSP1 in these solvents is approximately 1, 25, and 30 mg/ml, respectively.

iFSP1 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, iFSP1 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. iFSP1 has a solubility of approximately 0.1 mg/ml in a 1:7 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

iFSP1 is an inhibitor of apoptosis inducing factor mitochondria-associated 2/ferroptosis suppressor protein 1 (AIFM2/FSP1).¹ It decreases cell viability in glutathione peroxidase 4 (GPX4) knockout, but not wild-type, Pfa1 and HT1080 cells overexpressing AIFM2/FSP1, an effect that can be blocked by the ferroptosis inhibitor liproxstatin-1 (Item No. 17730). iFSP1 (3 μM) sensitizes a variety of human cancer cell lines to the ferroptosis inducer (1S,3R)-RSL3 (Item No. 19288).

Reference

1. Doll, S., Freitas, F.P., Shah, R., *et al.* FSP1 is a glutathione-independent ferroptosis suppressor. *Nature* (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, 03/20/2024

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