

PRODUCT INFORMATION



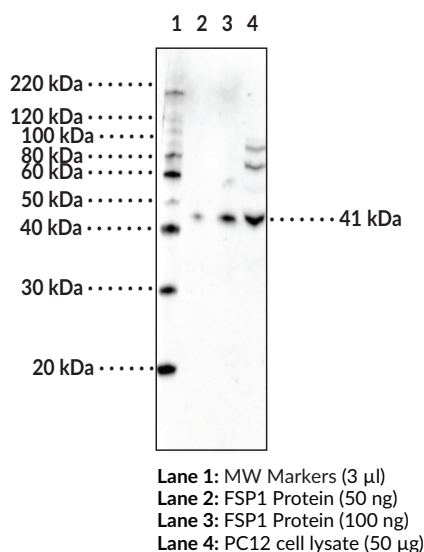
FSP1 Polyclonal Antibody

Item No. 31406

Overview and Properties

| | |
|----------------------------|---|
| Contents: | This vial contains 500 µg of protein A-affinity purified polyclonal antibody. |
| Synonyms: | AIFM2, Apoptosis-inducing Factor Homologous Mitochondrion-associated Inducer of Death, Apoptosis-inducing Factor Mitochondria-associated 2, Ferroptosis Suppressor Protein 1, p53-Responsive Gene 3 Protein, PRG3 |
| Immunogen: | Full-length human recombinant FSP1 |
| Cross Reactivity: | (+) FSP1 |
| Species Reactivity: | (+) Human, mouse |
| Uniprot No.: | Q9BRQ8 |
| Form: | Liquid |
| Storage: | -20°C (as supplied) |
| Stability: | ≥3 years |
| Storage Buffer: | PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide |
| Host: | Rabbit |
| Applications: | ELISA and Western blot (WB); the recommended starting dilution for ELISA and WB is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically. |

Image



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, 12/13/2023

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

PRODUCT INFORMATION



Description

Ferroptosis suppressor protein 1 (FSP1), also known as apoptosis-inducing factor mitochondria-associated 2 (AIFM2), is a flavoprotein and NAD(P)H-dependent oxidoreductase that inhibits ferroptosis in a glutathione peroxidase 4 (GPX4) and glutathione-independent manner.¹⁻³ It is encoded by *AIFM2* in humans and is composed of a short N-terminal hydrophobic region followed by a flavin adenine dinucleotide-dependent oxidoreductase domain.² FSP1 localizes to the plasma membrane in a myristoylation-dependent manner, where it reduces coenzyme Q₁₀ (CoQ₁₀; Item No. 11506) to CoQ₁₀H₂. Reduced CoQ₁₀ functions as a radical-trapping antioxidant and inhibits lipid peroxidation. *AIFM2* expression positively correlates with resistance to GPX4 inhibitors, including (1S,3R)-RSL3 (Item No. 19288), ML-210 (Item No. 23282), and ML-162 (Item No. 20455), in cancer cell lines. Withdrawal of the ferroptosis inhibitor ferrostatin-1 (Item No. 17729) reduces tumor growth in an *Fsp1/Gpx4* double knockout, but not a *GPX4* single knockout, H460 lung cancer mouse xenograft model, indicating that FSP1 confers resistance to ferroptotic cell death. Cayman's FSP1 Polyclonal Antibody can be used for ELISA and Western blot (WB) applications. The antibody recognizes FSP1 at 41 kDa from human and mouse samples.

References

1. Doll, S., Freitas, F.P., Shah, R., *et al.* FSP1 is a glutathione-independent ferroptosis suppressor. *Nature* **575(7784)**, 693-698 (2019).
2. Bersuker, K., Hendricks, J., Li, Z., *et al.* The CoQ oxidoreductase FSP1 acts parallel to GPX4 to inhibit ferroptosis. *Nature* **575(7784)**, 688-692 (2019).
3. Marshall, K.R., Gong, M., Wodke, L., *et al.* The human apoptosis-inducing protein AMID is an oxidoreductase with a modified flavin cofactor and DNA binding activity. *J. Biol. Chem.* **280(35)**, 30735-30740 (2005).