

FOLR1 (human) monoclonal antibody (548908)

The protein encoded by this gene is a member of the folate receptor (FOLR) family. Members of this gene family have a high affinity for folic acid, several reduced folic acid derivatives, and mediate delivery of 5-methyltetrahydrofolate to the interior of cells. Human FOLR1 shares 83% amino acid sequence identity with mouse and rat FOLR1. FOLR1 is up-regulated on many carcinomas, such as ovarian, and predominantly expressed on epithelial cells.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-ABS378-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable one month reconstituted under sterile conditions at 4°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Folate-binding protein, gp38, FRα, Folate Receptor α, Folate Receptor 1
Application	ELISA, Flow Cytometry, ICC, WB
Clone	548908
Crossreactivity	Does not cross-react with recombinant human FOLR2, 3 or 4 in direct ELISAs.
Formulation	Lyophilized from a 0.2 µm filtered solution in PBS with 5% Trehalose.
Host	Mouse
Immunogen	Chinese hamster ovary cell line CHO derived recombinant human FOLR1 Arg25Met233.
Isotype	IgG1
Purity Detail	Protein A/G purified from hybridoma culture supernatant.
Recommendation Dilutions/Conditions	ELISA (0.1-0.4 µg/ml (Detection), 2-8 µg/mL (Capture))Flow Cytometry (0.25 µg/106 cells)Immunocytochemistry (8-25 µg/mL)Western Blot (2 µg/mL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Reconstitution	Reconstitute at 0.5 mg/mL in sterile PBS.
Species Reactivity	Human
Technical Info / Product Notes	37 – 42 kDa protein
UniProt ID	P15328
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)