

FOLR2 recombinant monoclonal antibody (001)

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

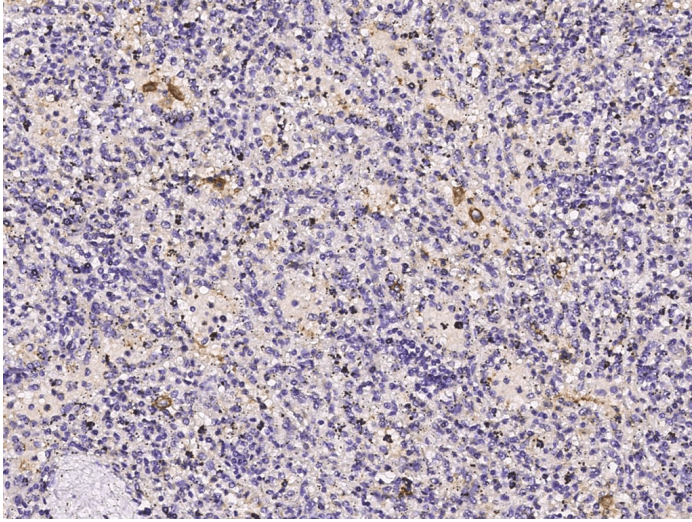
Ordering Information

[Order Online »](#)

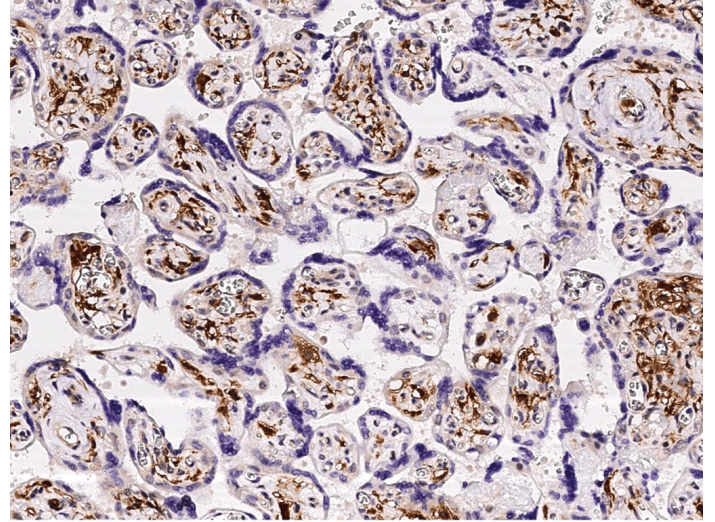
ENZ-ABS975-0100	100µl
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunochemical staining of human FOLR2 in human spleen with rabbit monoclonal antibody at 1:5000 dilution, formalin-fixed paraffin embedded sections.



Immunochemical staining of human FOLR2 in human placenta with rabbit monoclonal antibody at 1:5000 dilution, formalin-fixed paraffin embedded sections.

Handling & Storage

Handling Protected from light. Avoid freeze-thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Folate receptor 2, Folate receptor β , Placental folate-binding protein

Application IHC (PS)

Clone 001

Formulation Liquid. 0.2 μ m filtered solution in PBS. Preservative free.

Host Rabbit

Immunogen Recombinant human FOLR2 protein.

Isotype IgG

Purity Detail Protein A purified.

Recommendation Dilutions/Conditions Immunohistochemistry (paraffin sections) (1:2'000-1:10'000) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity Human

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

.



Last modified: September 17, 2024

ENZO LIFE SCIENCES

Phone: 800.942.0430
info-
usa@enzolifesciences.com

European Sales Office

ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-
eu@enzolifesciences.com

Belgium, The Netherlands

& Luxembourg
Phone: +32 3 466 0420
info-
be@enzolifesciences.com

France

Phone: +33 472 440 655
info-
fr@enzolifesciences.com

Germany

Phone: +49 7621 5500 526
info-
de@enzolifesciences.com

UK & Ireland

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