

FABP5 polyclonal antibody

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

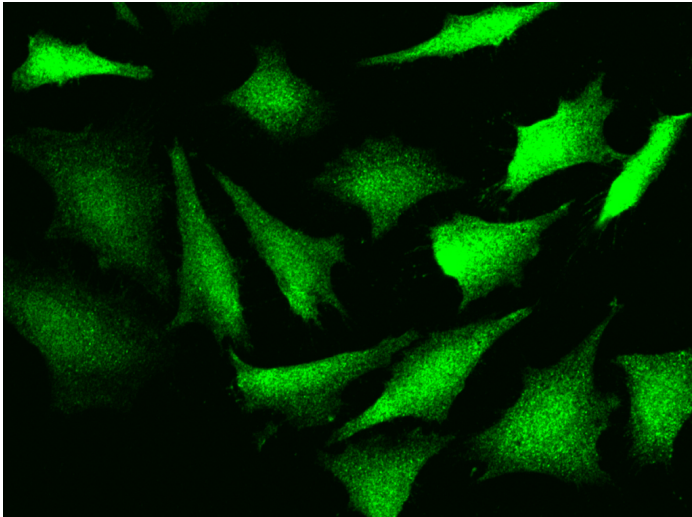
Ordering Information

[Order Online »](#)

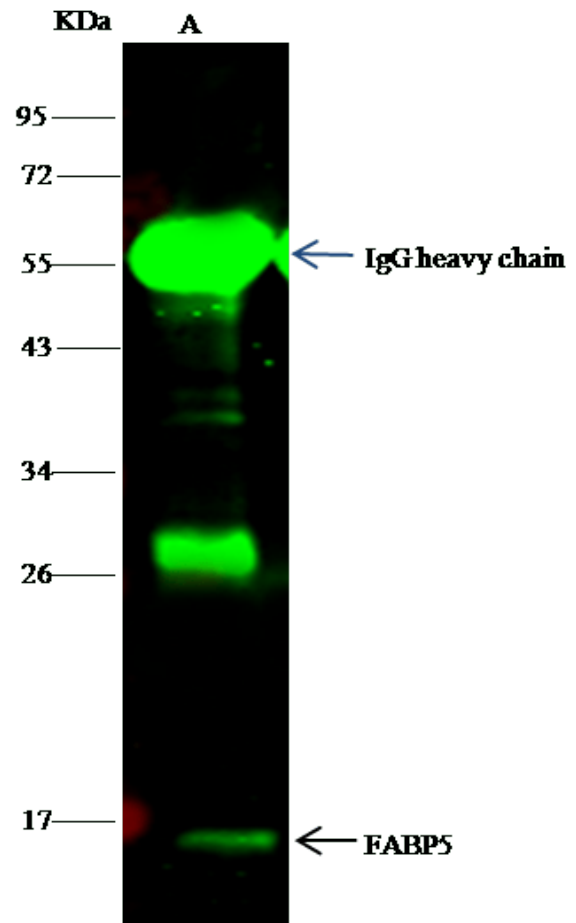
ENZ-ABS994-0100	100µl
-----------------	-------

Manuals, SDS & CofA

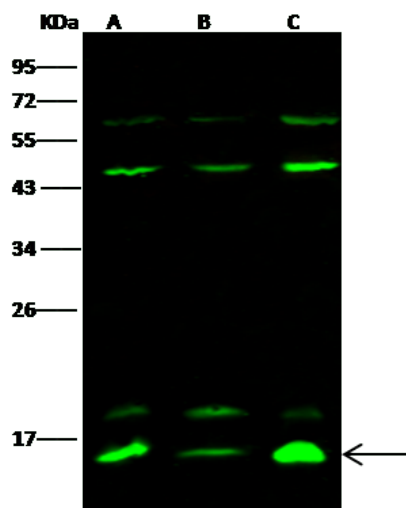
[View Online »](#)



Immunofluorescence staining of FABP5 in HeLa cells. Cells were fixed with 4% PFA, permeabilized with 0.3% Triton X-100 in PBS, blocked with 10% serum, and incubated with rabbit anti-human FABP5 polyclonal antibody (1:1000) at 4 °C overnight. Then cells were stained with the Alexa Fluor®488-conjugated Goat Anti-rabbit IgG secondary antibody (green). Positive staining was localized to cytoplasm and nucleus.



FABP5 was immunoprecipitated using: Lane A: 0.5 mg A431 Whole Cell Lysate, 4 µL anti-FABP5 rabbit polyclonal antibody and 15 µL of 50 % Protein G agarose. Primary antibody: Anti-FABP5 rabbit polyclonal antibody, at 1:100 dilution. Secondary antibody: DyLight 800-labeled antibody to rabbit IgG (H+L), at 1:5000 dilution. Developed using the Odyssey technique. Performed under reducing conditions. Observed band size: 15 kDa.



Anti-FABP5 rabbit polyclonal antibody at 1:500 dilution.
Lane A: HeLa Whole Cell Lysate Lane B: A431 Whole
Cell Lysate Lane C: HepG2 Whole Cell Lysate.
Lysates/proteins at 30 μ g per lane. Secondary Goat
Anti- Rabbit IgG H&L (Dylight 800) at 1/10000 dilution.
Developed using the Odyssey technique. Performed
under reducing conditions. Observed band size:15 kDa.

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	Fatty acid-binding protein 5, E-FABP, Epidermal-type fatty acid-binding protein
Application	ELISA, ICC, IF, IP, WB
Formulation	Liquid. 0.2 µm filtered solution in PBS. Preservative free.
Host	Rabbit
Immunogen	Recombinant human FABP5/E-FABP protein.
Isotype	IgG
Purity Detail	Protein A and antigen affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:500-1:1'000)ELISA (1:5'000-1:10'000)Immunocytochemistry (1:300-1:10'000)Immunofluorescence (1:300-1:10'000)Immunoprecipitation (4-8 µl/ml of lysate)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](#)



ENZO LIFE SCIENCES,
INC.
September 17, 2024
info-
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(EU) 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