

α -Fetoprotein recombinant monoclonal antibody (003)

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

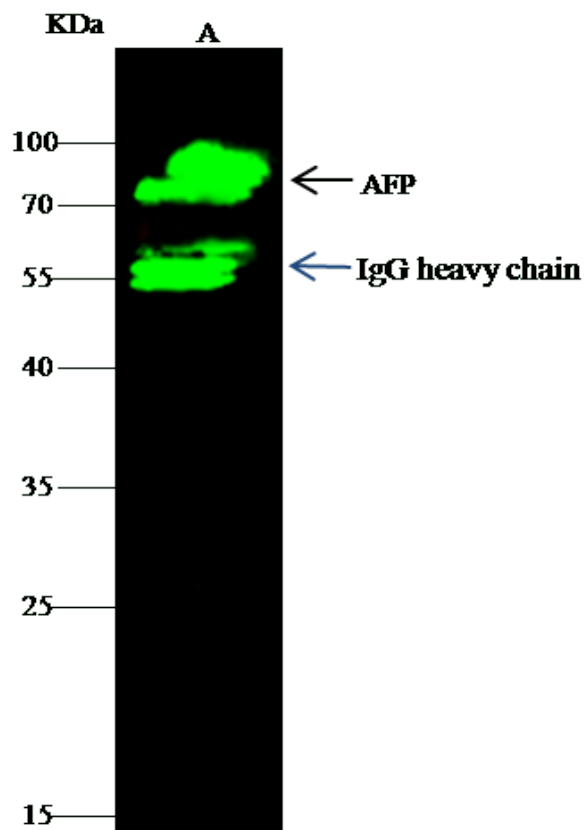
Ordering Information

[Order Online »](#)

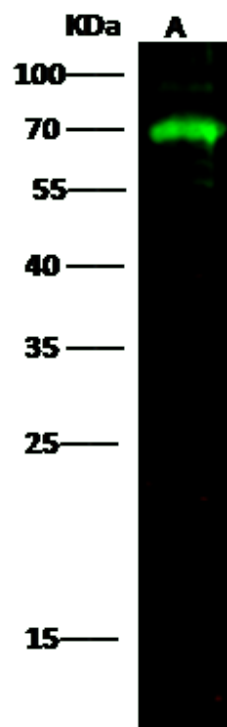
ENZ-ABS990-0100	100 μ l
-----------------	-------------

Manuals, SDS & CofA

[View Online »](#)



AFP was immunoprecipitated using: Lane A: 0.5 mg HepG2 Whole Cell Lysate, 2 μ L anti-AFP rabbit monoclonal antibody and 60 μ g of Immunomagnetic beads Protein G. Primary antibody: Anti-AFP rabbit monoclonal 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: 68 kDa.



Anti-AFP rabbit monoclonal antibody at 1:500 dilution. Lane A: HepG2 Whole Cell Lysate. Lysates/proteins at 30 μ g per lane. Secondary Goat Anti-Rabbit IgG H&L (Dylight800) at 1/10000 dilution. Developed using the Odyssey technique. Performed under reducing conditions. Observed band size: 69 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 AFP

Application ELISA, IP, WB

Clone 003

Formulation Liquid. 0.2 µm filtered solution in PBS. Preservative free.

Host Rabbit

Immunogen Recombinant human AFP / alpha-fetoprotein. protein.

Isotype IgG

Purity Detail Protein A purified.

Recommendation ELISA (1:5'000-1:10'000)Western Blot (1:500-1:2'000)Immunoprecipitation (0.5-2 µl/mg
Dilutions/Conditions 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.
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-
fr@enzolifesciences.com](mailto:info-fr@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)