

Angiopoietin-2 recombinant monoclonal antibody (117)

This monoclonal antibody detects Angiopoietin-2 (ANG2), a key regulator of angiogenesis and vascular remodeling.

ANG2 is a ligand of the Tie2 receptor and plays a critical role in blood vessel formation and remodeling. It is involved in vascular permeability and is frequently studied in cancer, inflammation, and cardiovascular diseases.

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

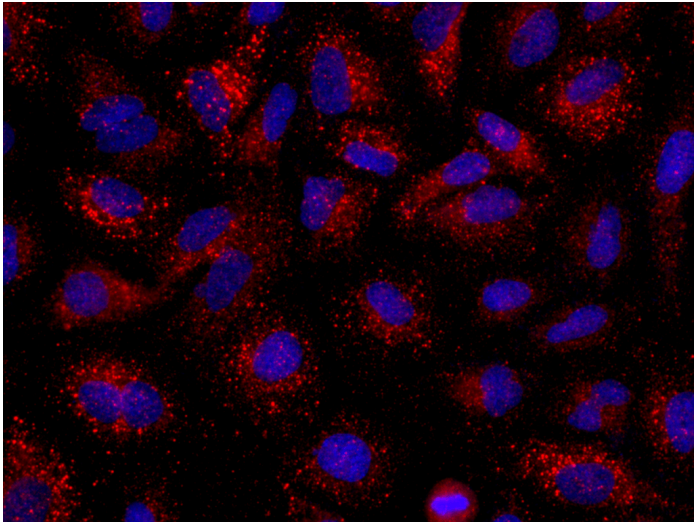
Ordering Information

[Order Online »](#)

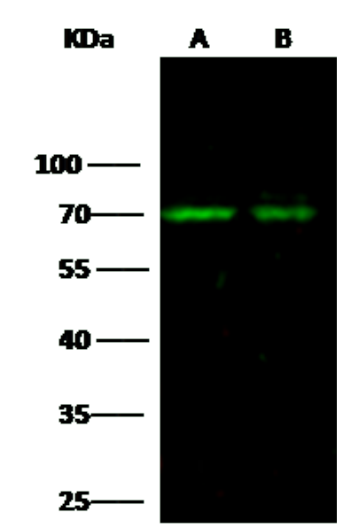
ENZ-ABS1066-0100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunofluorescence staining of Human Angiopoietin-2 in A549 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 Ang2 monoclonal antibody (1:300) at 4°C overnight. Then cells were stained with the Alexa Fluor® 549-conjugated Goat Anti-rabbit IgG secondary antibody (red) and counterstained with DAPI (blue).



Western blot analysis of various lysates using Angiopoietin-2 rabbit monoclonal antibody (ENZ-ABS1066) at 1:500 dilution. Secondary antibody: Goat Anti-Rabbit IgG H&L (Dylight800) at 1/10000 dilution. Lane A: HepG2 Whole Cell Lysate, Lane B: A549 Whole Cell Lysate. Lysates/proteins at 30 µg per lane.

Handling & Storage

Handling Avoid freeze-thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Ang2, ANGPT2

Application ICC, IF, IP, WB

Clone 117

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

Host Rabbit

Immunogen Recombinant human Angiopoietin-2.

Isotype IgG

Purity Detail Protein A purified.

Recommendation Immunocytochemistry/Immuofluorescence (1:100-1:500)

Dilutions/Conditions Western Blot (1:500-1:1'000)

Immunoprecipitation (0.2-1µL/mg 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-
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)