

IHC universal negative control reagent

IHC universal negative control reagent is specifically designed to work with rabbit, mouse and goat antibodies. The reagent has been titrated to work with polymer-based secondary systems in manual and automated IHC platforms.

Negative controls are used to determine the specificity of an antibody. No staining should be observed when the primary antibody is omitted. Minimal non-specific staining should be detected when using the IHC universal negative control reagent.

Citations: 12 [View Online »](#)

Ordering Information [Order Online »](#)

ADI-950-231-0025	25ml
------------------	------

Manuals, SDS & CofA [View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Formulation

Liquid. Negative control reagent- mixture of purified goat, mouse and rabbit immunoglobulins (IgGs).

Technical Info / Product Notes

Interpretation of the results will be the sole responsibility of the user. Source is from purified goat, mouse and rabbit immunoglobulins (IgGs) and total IgG concentration is 1.5 ug/ml.

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)