

Goat anti-rat IgG, polyclonal antibody (HRP conjugate)

Specificity: Based on immunoelectrophoresis and/or ELISA, the antibody reacts with whole molecule rat IgG. It also reacts with the light chains of other rat immunoglobulins. No antibody was detected against non-immunoglobulin serum proteins. The antibody has been tested by ELISA and/or solid-phase absorbed to ensure minimal cross-reaction with human, mouse, bovine, horse and rabbit serum proteins, but it may cross-react with immunoglobulins from other species.

Immunoglobulin G (IgG) is an immunoglobulin monomer consisting of two (gamma) heavy chains and two light chains. Each IgG molecule contains two antigen binding domains and a single effector (Fc) domain.

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

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-SAB-220-0500	0.5ml
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Warning that sodium azide as a preservative will substantially inhibit the enzyme activity of HRP.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	ELISA, ICC, IHC, WB
Formulation	Liquid. In 0.005M sodium phosphate, pH 7.6, containing 0.125M sodium chloride, 7.5mg/ml BSA (IgG-free, protease-free), and 50% glycerol.
Host	Goat
Immunogen	Purified rat IgG (whole molecule).
Purity Detail	Immunoaffinity purified.
Quality Control	The antibody has been tested by ELISA and/or solid-phase absorbed to ensure minimal cross-reaction with human, mouse, bovine, horse and rabbit serum proteins, but it may cross-react with immunoglobulins from other species.
Recommendation Dilutions/Conditions	ELISA (1:5,000-1:100,000)Immunocytochemistry (1:500-1:5,000)Immunohistochemistry(1:500-1:5,000)Western Blot (1:10,000-1:200,000, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Rat

Specificity

Based on immunoelectrophoresis and/or ELISA, the antibody reacts with whole molecule rat IgG. It also reacts with the light chains of other rat immunoglobulins. No antibody was detected against non-immunoglobulin serum proteins.

Worry-free Guarantee

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

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)