

Goat anti-rabbit IgG, polyclonal antibody (AP conjugate)

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. Assay Designs offers isotyping kits and antibodies in a variety of formats for the detection of human, mouse, rabbit, rat, sheep, and goat IgG.

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

Citations: 10

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-SAB-301-J	1ml
---------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ELISA, WB

Formulation Liquid. In 0.005M TRIS-hydrogen chloride, pH 8.0, containing 0.125M sodium chloride, 7.5mg/ml BSA (IgG-free, protease-free), 50% glycerol, and 0.025M sodium azide.

Host Goat

Immunogen Purified rabbit IgG (whole molecule).

Purity Detail Immunoaffinity purified.

Quality Control Goat antiserum immunoaffinity purified to isolate specific anti-rabbit IgG antibodies then adsorbed with human IgG and conjugated to alkaline phosphatase.

Recommendation
Dilutions/Conditions ELISA (1:2,500-25,000)Western Blot (1:2500-25,000, colorimetric)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Rabbit

Specificity Based on IEP and/or ELISA, reacts with whole molecule rabbit IgG, and also the light chains of other rabbit immunoglobulins. No antibody detected against non-immunoglobulin serum proteins. Antibody tested by ELISA and/or solid-phase adsorbed to ensure minimal cross-reaction with human serum proteins.

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)