

Goat anti-mouse IgG F(ab')₂, polyclonal antibody (HRP 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: 145

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-SAB-100-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, 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 Fab fragment of mouse IgG.

Purity Detail Immunoaffinity purified.

Quality Control The antibody has been tested by ELISA and/or solid-phase adsorbed to ensure minimal cross-reaction with human, bovine and horse serum proteins.

Recommendation ELISA (1:2,500-1:50,000)Immunohistochemistry (1:250-1:2500)Western Blot (1:2500-1:50,000 colorimetric, 1:5,000-1:100,000 ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Mouse

Specificity Based on immunoelectrophoresis and/or ELISA, the antibody reacts with the F(ab')₂/Fab portion of mouse IgG, and also with the light chains of other mouse immunoglobulins. No antibody was detected against the Fc portion of mouse IgG or against other non-immunoglobulin serum proteins.

Technical Info / Product Notes **Cited samples:**
[For an overview on cited samples please click here.](#)

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)