

# Goat anti-mouse IgG F(ab')<sub>2</sub>, 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: 149

[View Online »](#)

## Ordering Information

[Order Online »](#)

ADI-SAB-100-J	1ml
---------------	-----

## Manuals, SDS & CofA

[View Online »](#)

**Long Term Storage**      -20°C

**Regulatory Status** RUO - Research Use Only

<b>Application</b>	ELISA, IHC, WB
--------------------	----------------

Host	Goat
------	------

<b>Immunogen</b>	Fab fragment of mouse IgG.
------------------	----------------------------

<b>Purity Detail</b>	Immunoaffinity purified.
----------------------	--------------------------

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

<b>Recommendation Dilutions/Conditions</b>	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

<b>Specificity</b>	Based on immunoelectrophoresis and/or ELISA, the antibody reacts with the F(ab') <sub>2</sub> /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](#)

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)