

Goat anti-mouse IgG, (HRP conjugate)

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

Citations: 49

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SA204-0100	100µg
----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. For long term storage, aliquot product into individual tubes and freeze at -20 or -70°C. Defrost samples as needed and store remainder at 4°C for stability up to several days. Avoid repeated freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	ELISA, WB
Crossreactivity	Does not cross-react with human, horse, goat, sheep, bovine, guinea pig, hamster, rat, rabbit, or chicken proteins.
Formulation	Liquid. In 20 mM potassium phosphate, pH 7.2, containing 150mM sodium chloride, 0.01% gentamicin, and 1% BSA as stabilizer.
Host	Goat
Immunogen	Mouse IgG.
Purity Detail	Affinity purified.
Recommendation Dilutions/Conditions	Western Blot (secondary Ab detection, 1:5,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Mouse
Specificity	Recognizes mouse IgG heavy and light chains (whole molecule).

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)