

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

Immunoglobulin M (IgM) is an immunoglobulin expressed initially on the surface of B-cells as a monomer of two mu heavy chains and two light chains. Following B-cell activation, IgM is secreted in its pentameric form, consisting of five IgM molecules joined at the base of the Fc domain by a J chain. IgM is the first Immunoglobulin expressed following infection.

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

Ordering Information

[Order Online »](#)

ADI-SAB-210-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 PBS, pH 7.6, containing 7.5mg/mL BSA and 50% glycerol.
Host	Goat
Immunogen	Purified rat IgM (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, bovine, and horse serum proteins, but it may cross-react with IgM from other species.
Recommendation Dilutions/Conditions	ELISA (1:2,500-50,000)Western Blot (1:5,000-100,000, 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 the heavy chain of rat IgM but not with rat IgG or the light chains of rat immunoglobulins. No antibody was detected against non-immunoglobulin 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)

Last modified: May 29, 2024