

# Goat anti-mouse 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](#).

Citations: 27

[View Online »](#)

## Ordering Information

[Order Online »](#)

ADI-SAB-110-J	1ml
---------------	-----

## Manuals, SDS & CofA

[View Online »](#)





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)