

IgM (mouse), ELISA kit

The IgM (mouse), ELISA kit is a colorimetric immunometric enzyme immunoassay kit with results in 90 minutes.

The IgM (mouse), EIA kit is a colorimetric immunometric enzyme immunoassay kit with results in 90 minutes. Absorbance is read at 450 nm. Ideal for determining the isotype and concentration of monoclonal antibodies.

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.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

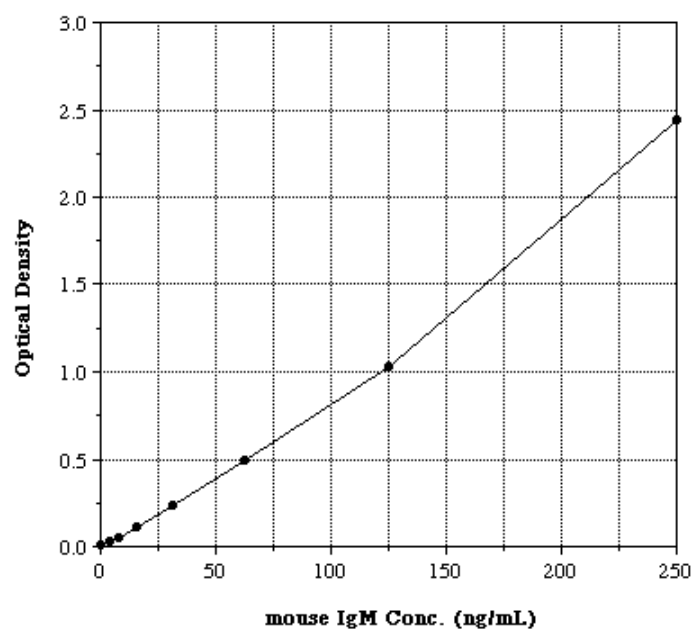
ADI-900-120

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Time savings - determine the isotype and concentration of monoclonal antibodies
- Quantitative - suitable in determining concentration
- Specific - no reactivity with other immunoglobulins
- Fast - results from up to 40 samples in duplicate in just 90 minutes



Handling & Storage

Use/Stability Store all components at 4°C, except standard at -20°C.

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of IgM in ascites fluid, culture supernatants, and serum of mouse origin.

Assay Time 90 minutes

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

Contents GxM IgG Microtiter plate, mouse IgM conjugate, Assay buffer 13 concentrate, Wash buffer concentrate, mouse IgGM standard, TMB Substrate, Stop solution 2

Crossreactivity Mouse IgM (100%), rat IgM (60%), human IgM (<0.4%), mouse IgG1 (<0.2%) and <0.10%: mouse IgG2a, mouse IgG2b, mouse IgG3

Sensitivity 0.60 ng/ml (range 3.91 – 250 ng/ml)

Species Reactivity Mouse

Wavelength 450 nm



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)