

IgG₁ (mouse), ELISA kit

Highly specific quantitative ELISA kit with results in 90min

The IgG₁ (mouse), EIA kit is a colorimetric immunometric enzyme immunoassay kit with results in 90 minute. Absorbance is read at 450 nm. Ideal for determining the isotype and concentration of monoclonal antibodies.

Immunoglobulin G (IgG) is an immunoglobulin monomer consisting of two (γ) heavy chains and two light chains. Each IgG molecule contains two antigen binding domains and a single effector (Fc) domain. We offer isotyping kits and antibodies in a variety of formats for the detection of human, mouse, rabbit, rat, sheep, and goat IgG.

Ordering Information

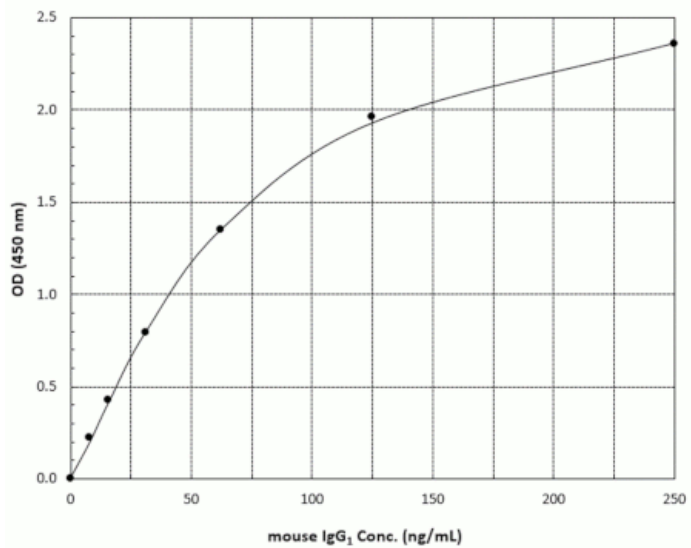
[Order Online »](#)

ADI-900-109A	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
- Results from up to 40 samples in duplicate in just 90 minutes



Handling & Storage

Use/Stability Store all components at 4°C, except the conjugate concentrate at -20°C.

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of IgG₁ in 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 IgG1 conjugate concentrate, Assay buffer 13 concentrate, Wash buffer concentrate, mouse IgG1 conjugate diluent, mouse IgG1 standard, TMB Substrate, Stop solution 2

Crossreactivity Mouse IgG1 (100%), rat IgG1 (0.90%), mouse IgG2b (0.21%) and <0.01%: human IgG1, mouse IgG2a, mouse IgG3, mouse IgM

Sensitivity 0.116 ng/mL (range 7.81 – 250 ng/ml)

Species Reactivity Mouse



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)