

IgG_{2b} (mouse), ELISA kit

The IgG_{2b} (mouse), ELISA kit is a colorimetric immunometric enzyme immunoassay kit with results in 90 minutes.

The IgG_{2b} (mouse), EIA kit is a colorimetric immunometric enzyme immunoassay kit with results in 90 minutes. 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

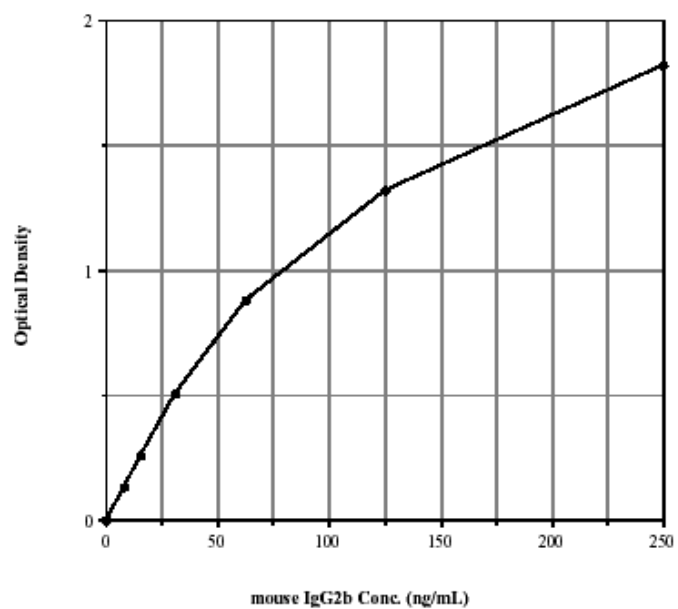
[Order Online »](#)

ADI-900-110	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 min



Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of IgG_{2b} 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 IgG2b conjugate, Assay buffer 13 concentrate, Wash buffer concentrate, mouse IgG2b standard, TMB Substrate, Stop solution 2

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

Sensitivity 1.05 ng/ml (range 7.81 – 250 ng/ml)

Species Reactivity Mouse

Wavelength 450 nm



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com