

# Mouse IgG Isotyping Kit ab273149

## Overview

**Product name** Mouse IgG Isotyping Kit

**Product overview** The Mouse IgG Isotyping Kit (ab273149) provides a rapid and easy method (one step ELISA) to characterize mouse monoclonal antibody isotypes in cell culture supernatants or purified antibodies preparations. The kit includes ready-to-use reagents necessary to analyze 12 samples in less than 30 minutes.

Anti-mouse antibodies specific to each of the common light and heavy chains are pre-coated in the wells. Samples are pipetted into microwells and Ig present in the sample are bound by the capture antibody. Then, a HRP (horseradish peroxidase) conjugated anti-mouse IgG (H+L) antibody is pipetted and incubated simultaneously with samples. After washing microwells in order to remove any non specific binding, the ready to use substrate solution (TMB) is added to microwells and color develops if the specific immunoglobulin is present in the sample. Color development is then stopped by addition of stop solution. Absorbance is measured at 450 nm.

## Properties

**Storage instructions** Store at +4°C. Please refer to protocols.

Components	1 x 96 tests	5 x 96 tests	1 x 96 tests
Detection antibody	1 x 12ml	1 x 60ml	1 x 12ml
Pre-coated microwells strips	1 x 96 tests	5 x 96 tests	1 x 96 tests
Sample Diluent	1 x 30ml	1 x 150ml	1 x 30ml
Stop solution	1 x 12ml	1 x 60ml	1 x 12ml
TMB Substrate	1 x 12ml	1 x 60ml	1 x 12ml

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

**Our Abpromise to you: Quality guaranteed and expert technical support**

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

#### **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors