

# Pro-Detect™ Rapid Antibody Isotyping Assay Kit - Mouse

Catalog Number A38550

Pub. No. MAN0017621 Rev. B



**WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from [thermofisher.com/support](https://www.thermofisher.com/support).

## Product description

The Thermo Scientific™ Pro-Detect™ Rapid Antibody Isotyping Assay Kit - Mouse (Cat. No. [A38550](#)) is a 5-minute lateral-flow assay offering ELISA-like sensitivity for the determination of mouse antibody classes and subclasses. To perform the assay, simply dip the Pro-Detect™ Rapid Antibody Isotyping Strips into a properly diluted tissue culture supernatant or mouse ascites sample.

Gold conjugates embedded in the membrane form specific, soluble complexes with the antibodies in the sample. These complexes migrate along the membrane and are captured by anti-isotype and class-specific antibodies impregnated in the membrane.

Results are displayed above the printed area on the strip as a red band, indicating the antibody isotype or subclass. A red band at the control (C) band region will always appear, confirming the test is functioning correctly.

## Contents and storage

Item	Storage
Pro-Detect™ Rapid Antibody Isotyping Strip 1 - Mouse	Store at 4°C. Do not freeze. After opening, store unused strips in the enclosed container containing desiccant.
Pro-Detect™ Rapid Antibody Isotyping Strip 2 - Mouse	
Pro-Detect™ Rapid Isotyping Dilution Buffer, 4 mL	

**Note:** Microtiter plates and test tubes are required to use the assay.

## Diute samples

**Note:** Proper sample dilution is essential for optimal results. If sample concentration is known, dilute to 100 ng/mL. For samples with unknown concentrations, follow the dilution instructions below.

1. Add 200 µL of Pro-Detect™ Rapid Isotyping Dilution Buffer to each culture tube.
2. Dilute 1:100 by adding 2 µL of fresh sample to the labeled culture tubes, then vortex to mix. For fluid samples containing antibodies at <10 µg/mL, dilute the sample 1:10 (instead of 1:100) by adding 22 µL of fluid sample.

## Perform the assay

1. Remove the container of lateral flow strips from refrigeration and fully equilibrate to room temperature.
2. Open the white tubes containing each of the lateral flow isotyping strips (Strip 1 and Strip 2) and add 150-200 µL of diluted sample to a well of a microtiter plate or test tube for each strip used.
3. Wait 5-10 minutes for colored bands to appear and immediately evaluate the results.

A successful test will result in a red band at the control (C) band region. A red band that develops towards the top or slightly above the printing for the class, subclass, or light chain type indicates the presence of that class, subclass, or light chain type in the fluid specimen tested.

**Note:** Keep all the test strips stored and sealed in desiccated tubes when not in use.

## Troubleshooting

Observation	Possible cause	Recommended action
No bands detected after applying sample.	Sample did not contain antibody.	Use a different test (e.g., ELISA or Western blot) to verify sample contains antibody . Verify correct assay strip is being used.
	Lateral flow strip was not sufficiently submerged in sample.	Apply test strip fully into sample well and ensure enough volume is present in the well to fully cover the white application tip of the strip.
Multiple bands detected.	Sample was not sufficiently diluted.	Dilute sample an additional 1:10 to 1:100 and re-test.
	Sample contained antibody from multiple sources.	Dilute sample an additional 1:10 to 1:100 and re-test. The darkest band is the isotype of the most abundant antibody in the sample.
	Ascites used as a source of antibody may have also contained antibodies from the host system.	Dilute sample an additional 1:10 to 1:100 and re-test. The darkest band is the isotype of the most abundant antibody in the sample.
	Myeloma cell line used in hybridoma production may have also secreted immunoglobulins.	Dilute sample an additional 1:10 to 1:100 and re-test. The darkest band is the isotype of the most abundant antibody in the sample.
	The test was allowed to proceed longer than 10 minutes.	Evaluate results within 10 minutes of adding the sample to the cassette well.

## Limited product warranty

Life Technologies Corporation and its affiliates warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have questions, contact Life Technologies at [www.thermofisher.com/support](http://www.thermofisher.com/support).



Thermo Fisher Scientific | 3747 N. Meridian Road | Rockford, Illinois 61101 USA

For descriptions of symbols on product labels or product documents, go to [thermofisher.com/symbols-definition](http://thermofisher.com/symbols-definition).

Revision history: Pub. No. MAN0017621 B

Revision	Date	Description
B	12 December 2024	The volume of Pro-Detect™ Rapid Isotyping Dilution Buffer was changed from 15 mL to 4 mL in the contents and storage table. Version numbering was changed to new format and reset to B in conformance with internal document control procedures.
A.0	22 May 2018	New document for Pro-Detect™ Rapid Antibody Isotyping Assay Kit - Mouse.

The information in this guide is subject to change without notice.

**DISCLAIMER:** TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

**Important Licensing Information:** These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2018-2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.