

Pro-Detect™ Rapid Antibody Isotyping Assay Kit - Human

Catalog Numbers A38552

Pub. No. MAN0017623 Rev. C



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 - Human (Cat. No. [A38552](#)) is a 5-minute lateral-flow assay with ELISA sensitivity for the determination of human antibodies. To perform the assay, simply dip the Pro-Detect™ Rapid Antibody Isotyping Strips into a properly diluted tissue-culture supernatant 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 printing on the strip as a red band, indicating the antibody isotype or subclass. A red colored band at the control (C) band region will always appear, confirming the test is functioning properly.

Contents and storage

Component	Storage
Pro-Detect™ Rapid Antibody Isotyping Strip 1 - Human	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 - Human	
Pro-Detect™ Rapid Isotyping Dilution Buffer, 4 mL	

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

Prepare samples

Note: Proper sample dilution is essential for optimal results.

Dilute samples as indicated.

For...	Do this...
Samples with known concentrations	Dilute samples to a concentration of 1 µg/mL to 5 µg/mL with Pro-Detect™ Rapid Isotyping Dilution Buffer. For example, to dilute a 10 µg/mL sample to 1 µg/mL, do the following. <ol style="list-style-type: none"> Add 200 µL of Pro-Detect™ Rapid Isotyping Dilution Buffer to a labeled culture tube. Add 22 µL of sample to the tube, then vortex to mix.
Samples with unknown concentrations	Dilute samples 1:100 with Pro-Detect™ Rapid Isotyping Dilution Buffer. <ol style="list-style-type: none"> Add 200 µL of Pro-Detect™ Rapid Isotyping Dilution Buffer to a labeled culture tube. Dilute the sample 1:100 by adding 2 µL of sample to the tube, then vortex to mix. <p>Note: Depending on the results, if the 1:100 dilution does not give a sensitive result, add more than 2 µL of sample.</p>

Perform the test

Before you begin, remove the container of lateral flow isotyping strips from refrigeration and fully equilibrate to room temperature.

1. Open the white tubes containing the isotyping strips (Strip 1 and Strip 2), then remove the required number of strips.

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

2. For each strip, add 150 µL to 200 µL of diluted sample to a well of a microtiter plate or test tube.
3. Dip one strip into each sample-containing well or tube, wait 5–10 minutes for colored bands to appear, then immediately evaluate the results.

A successful test will result in a red band at the "C" or control location. A red line 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.

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.
	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. If you have questions, contact Life Technologies at 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.

Revision history: Pub. No. MAN0017623 C

Revision	Date	Description
C	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 C in conformance with internal document control procedures.
B.0	27 May 2021	Minimal corrections done in the document as requested by product manager and published.
A.0	11 June 2018	New document for Pro-Detect™ Rapid Antibody Isotyping Assay Kit - Human.

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.