

## Product datasheet

# TMRM Assay Kit (Mitochondrial Membrane Potential) ab228569

★★★★★ 1 Abreviews 1 Image

### Overview

<b>Product name</b>	TMRM Assay Kit (Mitochondrial Membrane Potential)
<b>Detection method</b>	Fluorescent
<b>Sample type</b>	Adherent cells, Suspension cells
<b>Assay type</b>	Cell-based
<b>Product overview</b>	TMRM Assay Kit (Mitochondrial Membrane Potential) ab228569 is designed to quantify changes in mitochondrial membrane potential in living cells using flow cytometry, fluorescence plate reader, or fluorescence microscopy. It uses the red-fluorescent probe TMRM (tetramethylrhodamine, methyl ester) that localizes in mitochondria and detects mitochondrial membrane depolarization.

#### Related assay kits

TMRE is the ethyl ester of tetramethylrhodamine and is a close equivalent of TMRM. TMRE is used about 3 times more than TMRM for mitochondrial membrane potential assays. TMRE-Mitochondrial Membrane Potential Assay Kit [ab113852](#) is our most popular assay kit for mitochondrial membrane potential analysis.

**Notes** TMRM staining is used to monitor mitochondrial function. When accumulated in a negatively charge polarized mitochondria, TMRM fluoresces orange. When mitochondrial membrane potential collapses in apoptotic or metabolically stressed cells, the TMRM reagent is dispersed throughout the cell cytosol and fluorescence levels drop dramatically.

This kit provides enough reagents to analyze 5 x 96 wells (500 samples).

**Platform** Microplate reader, Fluor. microscope, Flow cyt.

### Properties

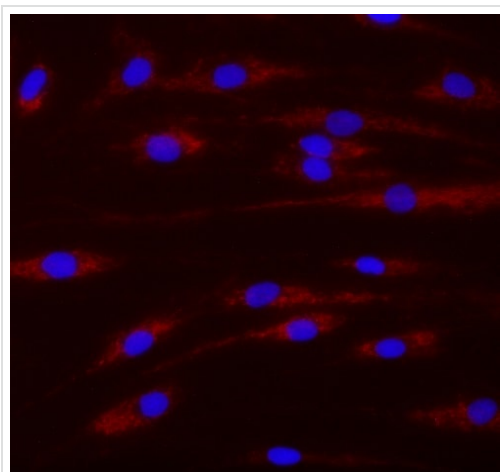
**Storage instructions** Please refer to protocols.

Components	1 kit
5X Live Cell Imaging Buffer	1 x 30ml
FCCP	1 x 25µl

Components	1 kit
TMRM Reagent	1 x 100µl

## Images

---



Example data

Human fibroblast cells stained using ab228569 TMRM Assay Kit (Mitochondrial Membrane Potential) (200 nM TMRM).

**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