

# Lipid Peroxidation (MDA) Assay Kit (Colorimetric) ab233471

[16 References](#) [2 Images](#)

### Overview

<b>Product name</b>	Lipid Peroxidation (MDA) Assay Kit (Colorimetric)
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Adherent cells, Suspension cells, Tissue Lysate
<b>Product overview</b>	Lipid Peroxidation (MDA) Assay Kit (Colorimetric) ab233471 enables researchers to detect MDA <u>without the heating steps</u> required by the TBARS assay conventionally used for MDA detection.

In this MDA assay, the MDA Color Reagent reacts with MDA to generate a blue color product which is measured at 695 nm with absorbance microplate readers. The assay is very fast and specific for MDA with little interference from other aldehydes.

Alternatively, see our popular [TBARS assay kit for MDA measurement ab118970](#).

MDA assay protocol summary for ab233471:

- add samples and standards to wells
- add MDA color reagent and incubate for 10-30 min at room temp
- add reaction solution and incubate for 30-60 min at room temp
- analyze with microplate reader

<b>Notes</b>	<p>Lipid peroxidation is characterized by the oxidative degradation of unsaturated fatty acids, phospholipids, glycolipids, cholesterol esters and cholesterol. Malondialdehyde (MDA) is one of the most commonly used biomarkers for lipid peroxidation.</p> <p>Running an MDA assay has historically relied on a reaction with thiobarbituric acid (the TBARS assay) to generate a product that can be measured colorimetrically at 532 nm or fluorimetrically at Ex/Em = 530/550 nm.</p> <p>However, the TBARS assay has quite a few limitations:</p> <ul style="list-style-type: none"> <li>- the reaction is not specific to MDA,</li> <li>- the TBA-MDA reaction needs be run under acidic conditions,</li> <li>- the TBARS assay needs be run under high temperature, commonly at 90-100 °C.</li> </ul>
--------------	--

<b>Tested applications</b>	<b>Suitable for:</b> Functional Studies
<b>Platform</b>	Microplate reader

## Properties

### Storage instructions

Store at -20°C. Please refer to protocols.

Components	200 tests
Dilution Buffer	1 x 10ml
MDA Color Reagent	1 vial
MDA Standard	1 vial
Reaction Solution	1 x 10ml

## Applications

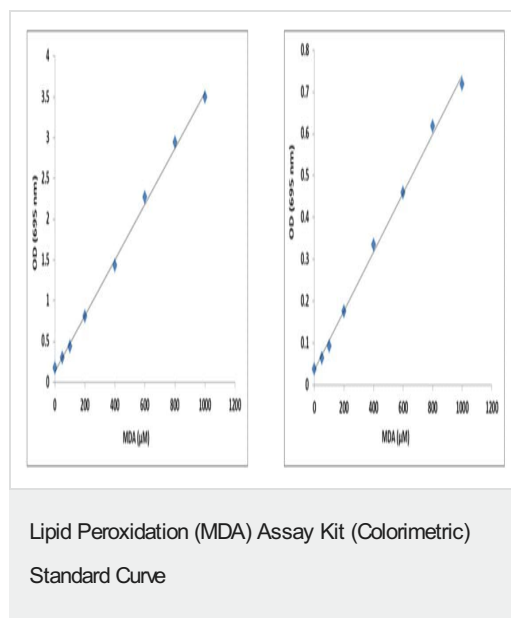
### The Abpromise guarantee

Our **Abpromise guarantee** covers the use of ab233471 in the following tested applications.

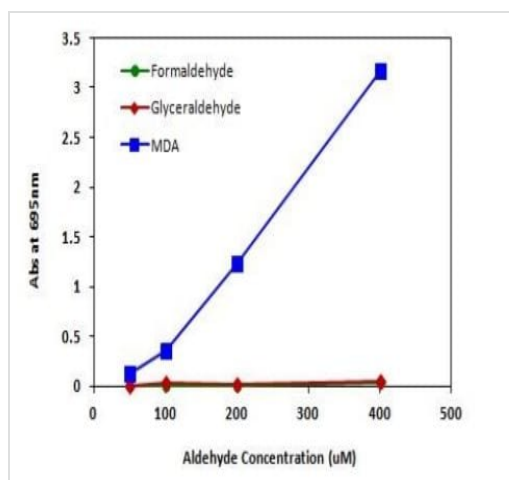
The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
Functional Studies		Use at an assay dependent concentration.

## Images



MDA dose response was measured with AB233471 on a 96-well clear bottom microplate using a SpectraMax microplate reader (Molecular Devices). (Pathcheck on (Left image); Pathcheck off (Right image))



Signal Comparison of MDA, Formaldehyde, and Glyceraldehyde

Signal Comparison of MDA, Formaldehyde, and Glyceraldehyde

**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