

# MDA Assay Kit (competitive ELISA) ab238537

## 24 References

### Overview

<b>Product name</b>	MDA Assay Kit (competitive ELISA)
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Serum, Plasma, Cell Lysate, Tissue Lysate
<b>Assay type</b>	Competitive
<b>Species reactivity</b>	<b>Reacts with:</b> Vertebrata
<b>Product overview</b>	MDA Assay Kit (competitive ELISA) (ab238537) is designed for the rapid detection and quantitation of Malondialdehyde (MDA)-protein adducts.

MDA is a widely accepted marker for oxidative stress. This kit will allow the quantitation of MDA adduct in protein samples determined by comparing its absorbance with that of a known MDA-BSA standard curve.

<b>Platform</b>	Microplate reader
-----------------	-------------------

### Properties

<b>Storage instructions</b>	Please refer to protocols.
-----------------------------	----------------------------

Components	96 tests
100X Conjugate Diluent	1 x 300µl
10X Wash Buffer	1 x 100ml
Anti-MDA Antibody (1000X)	1 x 10µl
Assay Diluent	1 x 50ml
MDA Conjugate	1 x 20µl
MDA-BSA Standard	1 x 50µl
Protein Binding Strip Well Plate	1 unit
Secondary Antibody, HRP Conjugate (1000X)	1 x 20µl

Components	96 tests
Stop Solution	1 x 12ml
Substrate Solution	1 x 12ml

**Relevance** Lipid peroxidation refers to the oxidative degradation of lipids and is a well-defined mechanism of cellular damage. The formation of lipid peroxidation products leads to spread of free radical reactions leading to cell damage.

**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