

## Product datasheet

# Aldehyde Assay Kit II (Colorimetric - Blue) ab219923

[1 Image](#)

### Overview

---

<b>Product name</b>	Aldehyde Assay Kit II (Colorimetric - Blue)
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Other biological fluids, Cell Lysate
<b>Assay type</b>	Quantitative
<b>Species reactivity</b>	<b>Reacts with:</b> Mammals, Other species
<b>Product overview</b>	Aldehyde Assay Kit II (Colorimetric - Blue) (ab219923) provides a sensitive mix-and-read method to detect aldehyde in a variety of samples such as cell lysates or biological fluids. This assay used a proprietary sensor that upon reacting with an aldehyde, generates a chromogenic product that can be detected by absorbance on a microplate reader at OD620-660 nm. This assay can detect as little as 3 $\mu$ M aldehyde (0.3 nmol/100 $\mu$ L assay volume).

The assay can be performed in a convenient 96-well or 384-well plate format and is easily adapted to automation without a separation step.

If your plate reader does not have the appropriate filter to detect the absorbance change, we recommend **Aldehyde Quantification Assay Kit (Colorimetric) (ab112113)**, which can be read at OD405 or OD550 nm.

**Notes** Very reactive aldehydes, namely 4-hydroxyalkenals, were first shown to be formed in autoxidizing chemical systems. It was subsequently shown that 4-hydroxyalkenals, particularly 4-hydroxynonenal, were formed in substantial amounts under biological conditions, i.e. during the peroxidation of lipids of liver microsomes incubated in the NADPH-Fe system. Many other aldehydes were also identified in peroxidizing liver microsomes or hepatocytes, e.g., alkanals, alk-2-enals, and 4-hydroxyalkenals.

Rapid and accurate measurement of aldehydes is an important task for biological and chemical research, as well as food industry and environmental pollution surveillance.

**Platform** Microplate reader

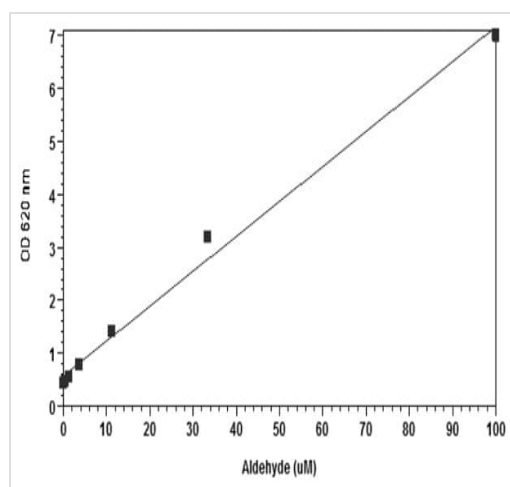
### Properties

---

**Storage instructions** Store at -20°C. Please refer to protocols.

Components	200 tests
Aldehyde Standard	1 vial
Assay Buffer	1 x 25ml
Blue Dye	2 units
Blue Enhancer	1 x 10ml

## Images



Aldehyde Quantification Assay Kit II (Colorimetric)  
(ab219923)

Aldehyde Quantification Assay Kit II (Colorimetric) (ab219923). Typical aldehyde standard calibration curve. Aldehyde dose response was measured in a white wall/clear bottom 96-well plate using a SpectraMax microplate reader. As low as 3  $\mu$ M aldehyde can be detected with 30 minutes incubation (n=3).

**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

