

Aldehyde Assay Kit (Fluorometric) ab138882

5 References 5 Images

Overview

Product name	Aldehyde Assay Kit (Fluorometric)
Detection method	Fluorescent
Sample type	Saliva, Urine, Plasma, Cell Lysate, Cell culture media
Assay type	Quantitative
Sensitivity	3 µM
Species reactivity	Reacts with: Mammals, Other species
Product overview	Abcam's Aldehyde Quantification Assay Kit (Fluorometric) (ab138882) is used for quantifying aldehydes at higher pH by using a proprietary fluorogenic dye that generates a strongly fluorescent product upon reacting with an aldehyde. This fluorimetric kit provides a sensitive mix-and-read method to detect as little as 0.3 nanomole of aldehyde/100 µL assay volume (3 µM). The assay can be performed in a convenient 96-well or 384-well microtiter-plate format and easily adapted to automation without a separation step. Its signal can be read by a fluorescence microplate reader at Ex/Em = 365/435 nm.
Notes	The formation, reactivity and toxicity of aldehydes originating from the peroxidation of lipids of cellular membranes have received great attention in recent years. Rapid and accurate measurement of aldehydes is an important task for biological research, chemical research, food industry and environmental pollution surveillance. There are a few reagents or assay kits available for quantifying the number of aldehydes. Most of the existing aldehyde test methods are based on separations either by the tedious and expensive HPLC-MS or GC-MS.
Platform	Microplate reader

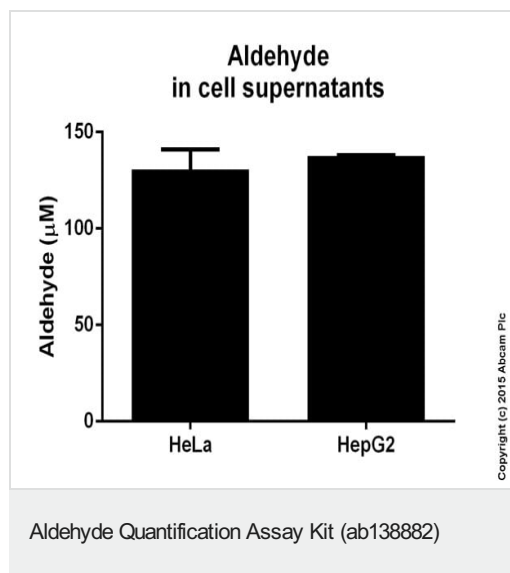
Properties

Storage instructionsStore at -20°C. Please refer to protocols.

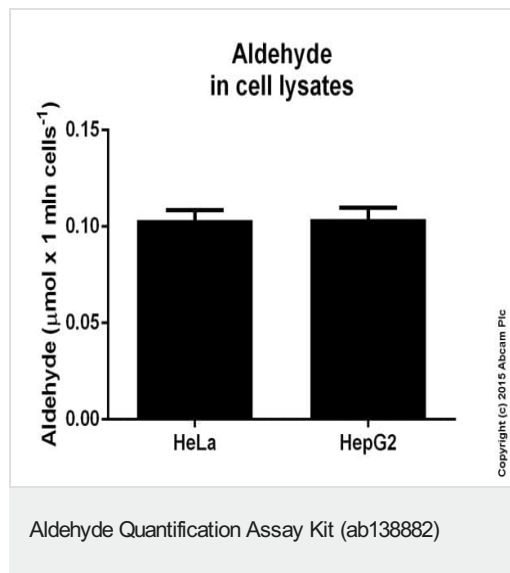
Components	200 tests
Aldehyde Blue Indicator	1 vial
Aldehyde Reaction Buffer	1 x 6ml
Aldehyde Standard	1 vial

Components	200 tests
Assay Buffer	1 x 30ml
DMSO	1 x 100µl

Images

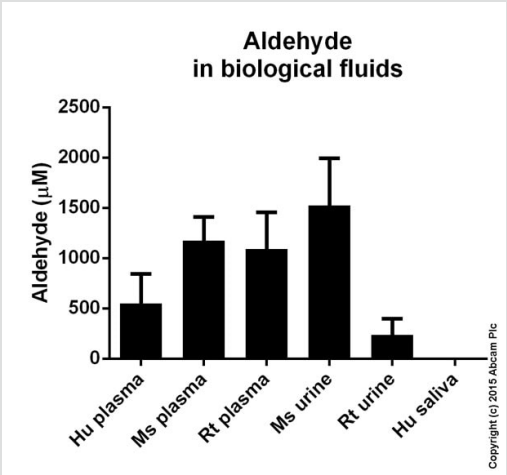


Aldehyde levels measured in cell supernatants (µM). Samples were diluted 2 fold.



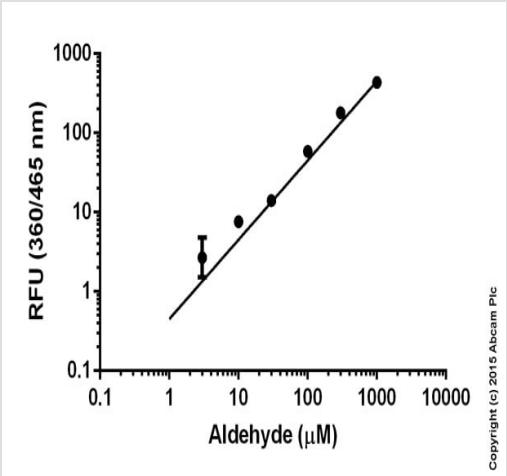
Aldehyde levels measured in cell lysates showing quantity (µmol) per 1 mln cells.

Samples with the concentration of 1e8 cells/mL (HepG2) and 1.28e8 cells/mL (HeLa) were used. Samples were diluted 100 fold.



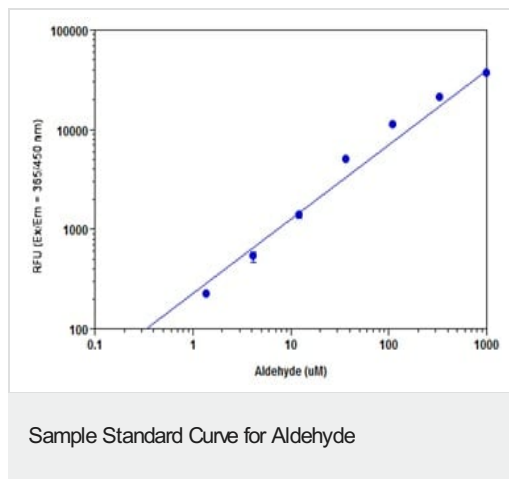
Aldehyde Quantification Assay Kit (ab138882)

Aldehyde levels measured in biological fluids (µM). Samples were diluted 1-10 fold.



Aldehyde Quantification Assay Kit (ab138882)

Standard curve: mean of duplicates (+/- SD) with background reads subtracted



Aldehyde dose response was measured in a solid black 96-well plate with ab138882 using a fluorescence microplate reader. As low as 3 μM of aldehyde can be detected with 15 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