

## AB308327 – Acetaldehyde Assay Kit (Fluorometric)

Determination of acetaldehyde in various sample types.

For research use only - not intended for diagnostic use.

PLEASE NOTE: With the acquisition of BioVision by Abcam, we have made some changes to component names and packaging to better align with our global standards as we work towards environmental-friendly and efficient growth. You are receiving the same high-quality products as always, with no changes to specifications or protocols.

For overview, typical data and additional information please visit: [Abcam Website](#)

### Storage and Stability

Store the kit at -20 °C, protected from light. Briefly centrifuge all small vials prior to opening. Read the entire protocol before performing the assay.

### Materials Supplied

Item	Quantity	Storage Condition
Assay Buffer XXVIII/ACH Assay Buffer	25 ml	RT
Recombinant ALDH2/ALDH2	50 µl	-20°C
Acetaldehyde/ACH Standard (0.1 M)	0.5 ml	RT
ALDH2 Substrate Mix/ACH Substrate Mix	1 vial	-20 °C
PicoProbe I/ACH Probe (Avoid light)	0.4 ml	RT

### Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully utilize this assay:

- Dounce Tissue Homogenizer
- 96-well white plate with flat bottom
- Multi-well spectrophotometer (Fluorescent plate reader)

### Reagent Preparation

Assay Buffer XXVIII/ACH Assay Buffer, Acetaldehyde/ACH Standard (0.1 M) and ACH Probe: Warm to room temperature (RT) before use.

Recombinant ALDH2/ALDH2: Keep on ice until it thaws completely. Divide into aliquots and store at -20 °C. Avoid repeated freeze/thaw cycles.

ACH Substrate Mix: Reconstitute the vial in 220 µl Assay Buffer XXVIII/ACH Assay Buffer. Pipette up and down to dissolve completely. Keep on ice and store at -20 °C.

### Non-Enzymatic Cell Harvesting Protocol

#### Sample Preparation:

##### For liquid samples (e.g., saliva)

Transfer 0.5 ml of liquid sample into an Eppendorf tube. Centrifuge at 10,000 x g for 20 min at RT. Transfer the supernatant to a new Eppendorf tube. Dilute the sample using Assay Buffer XXVIII/ACH Assay Buffer, if necessary.

##### For solid or food samples (e.g., banana)

Weigh ~100 mg of sample and transfer to an Eppendorf tube. Add 0.5 ml of cold Assay Buffer XXVIII/ACH Assay Buffer into the tube and homogenize the sample using Dounce tissue homogenizer for 5 min. Centrifuge at 10,000 x g for 20 min at RT. Transfer the supernatant to a new Eppendorf tube. Dilute the sample using Assay Buffer XXVIII/ACH Assay Buffer, if necessary. Add 50 µl of the diluted supernatant into two parallel wells of a 96-well white plate labeled as Sample and Sample Background Control.

#### Δ Note:

1. For Unknown Samples, we recommend doing a pilot experiment and testing several doses to ensure that the readings are within the linear range of the Standard Curve.
2. Endogenous compounds may interfere with the reaction. To ensure accurate determination of ACH in the test samples, we recommend spiking samples with a known amount of Acetaldehyde/ACH Standard (50 µM).

### Acetaldehyde/ACH Standard Preparation:

Mix 10 µl of 0.1 M Acetaldehyde/ACH Standard with 990 µl Assay Buffer XXVIII/ACH Assay Buffer to generate 1 mM Acetaldehyde/ACH Standard. Further, mix 200 µl of 1 mM Acetaldehyde/ACH Standard with 800 µl of Assay Buffer XXVIII/ACH Assay Buffer to generate 200 µM Acetaldehyde/ACH Standard. Perform a 2-fold serial dilution of 200 µM Acetaldehyde/ACH Standard by mixing 0.5 ml of 200 µM Acetaldehyde/ACH Standard with 0.5 ml of Assay Buffer XXVIII/ACH Assay Buffer to generate 100 µM of Acetaldehyde/ACH Standard. Perform a 2-fold serial dilution of 100 µM Acetaldehyde/ACH Standard and 50 µM Acetaldehyde/ACH Standard to generate 50 µM and 25 µM of Acetaldehyde/ACH Standards respectively. Add 50 µl of 25, 50 and 100 µM of Acetaldehyde/ACH Standards into the desired wells of a 96-well white plate. Add 50 µl Assay Buffer XXVIII/ACH Assay Buffer to the 0 µM Acetaldehyde/ACH Standard well.

### Reaction Mix Preparation:

Mix enough reagents for the number of assays to be performed. For each well, prepare 50 µl Mix containing:

	Reaction Mix	Background Mix
Assay Buffer XXVIII/ACH Assay Buffer	45.5 µl	46 µl
Recombinant ALDH2/ALDH2	0.5 µl	---
ACH Substrate Mix	2 µl	2 µl
ACH Probe	2 µl	2 µl

Mix well and add 50 µl of Reaction Mix to all Standard and Sample wells and 50 µl of Background Mix to Sample Background Control wells respectively. Mix well.

### Measurement:

Measure the fluorescence at Ex/Em= 535/587 nm in kinetic mode at 25 °C for 30-60 min. The Acetaldehyde/ACH Standard Curve can be read in endpoint mode (at the end of incubation time).

### Calculation:

Subtract the 0 Standard reading from all Standard readings and plot the Acetaldehyde/ACH Standard Curve. For Samples and Sample Background Control wells, choose any two time points within the linear range of the curve (t1 & t2) and obtain the corresponding readings (RFU1 and RFU2). Calculate ΔRFUS for Samples, i.e. (ΔRFUS = RFUS2 – RFUS1) and ΔRFUSBC for

Sample Background Control ( $\Delta\text{RFUSBC} = \text{RFUSBC2} - \text{RFUSBC1}$ ). Then subtract the  $\Delta\text{RFUSBC}$  from  $\Delta\text{RFUS}$  to get the corrected Sample readings. Apply the corrected Sample readings to the Acetaldehyde/ACH Standard Curve to get C  $\mu\text{M}$  of ACH. Calculate the ACH concentration ( $\mu\text{g/g}$ ) in food samples as shown below

$$\text{ACH } (\mu\text{g/g}) = \text{C} \times \text{V} \times \text{MW} \times \text{D} / \text{W}$$

Where: **C** is the concentration of ACH from the Standard Curve (in  $\mu\text{M}$ )

**V** is the total sample volume (in L)

**MW** is the molecular weight of ACH (44 g/mol)

**D** is the sample dilution factor (if applicable, D = 1 for undiluted samples)

**W** is the weight of sample used (in g)

### Technical Support

Copyright © 2023 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

[www.abcam.com/contactus](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)