

## ab204695 – Adenosine Deaminase (ADA) Activity Assay Kit (Fluorometric)

For the detection of ADA activity.

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:

<http://www.abcam.com/ab204695>

### Storage and Stability

On receipt entire assay kit should be stored at -20°C, protected from light. Upon opening, use kit within 6 months.

### Materials Supplied

Item	Quantity	Storage Condition
10X ADA Assay Buffer/ADA Assay Buffer (10x)	25 mL	-20°C
ADA Converter	1 vial	-20°C
Converter Enzyme VIII/ADA Developer	1 vial	-20°C
ADA Positive Control	1 vial	-20°C
OxiRed Probe/ADA Probe	200 µL	-20°C
ADA Substrate	500 µL	-20°C
Inosine Standard/Inosine Standard (10 mM)	100 µL	-20°C

### Materials Required, Not Supplied

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

- 96-well plate with flat bottom. White plate is preferred for this assay.
- Fluorescence microplate reader
- Protease Inhibitor Cocktail
- Dounce homogenizer

### Reagent Preparation

- Before using the kit, spin the tubes prior to opening.

10X ADA Assay Buffer/ADA Assay Buffer (10x): Make 1x Assay Buffer by adding one part 10X ADA Assay Buffer/10x Assay Buffer to nine parts deionized water. Store at -20°C or 4°C. Bring to 37°C before use.

ADA Converter and Converter Enzyme VIII/ADA Developer: Reconstitute each with 210 µl ADA Assay Buffer and mix gently by pipetting. Briefly centrifuge to collect the contents at the bottom of the tube. Aliquot and store at -20°C. Avoid repeated freeze/thaw.

ADA Substrate: Aliquot and store at -20°C. Avoid repeated freeze/thaw.

ADA Positive Control: Reconstitute with 22 µl ADA Assay Buffer and mix gently by pipetting. Briefly centrifuge to collect the contents at the bottom of the tube. Aliquot and store at -20°C. Avoid repeated freeze/thaw.

### Assay Protocol

- Equilibrate all materials and prepared reagents to room temperature prior to use.

### Sample preparation:

1. Rinse tissue and transfer ~100 mg of fresh or frozen tissue (stored at -80°C) to a pre-chilled homogenizer.
2. Add 300 µl cold ADA Assay Buffer containing protease inhibitor cocktail (not provided) and thoroughly homogenize tissue on ice.
3. Transfer the tissue homogenate to a cold microfuge tube.
4. To prepare cell extract, add 150-300 µl cold ADA Assay Buffer containing protease inhibitor cocktail (not provided) to 1-5 x 10<sup>6</sup> fresh or frozen cells and pipette several times to disrupt the cells.
5. Transfer cell homogenate including cell debris to a cold microfuge tube and agitate on a rotary shaker at 4°C for at least 15 min.
6. Centrifuge the tissue or cell homogenate at 16,000 X g, 4°C for 10 min.
7. Transfer the clarified supernatant to a fresh pre-chilled tube & store on ice. Use lysates immediately to assay ADA activity.

**Δ Note:** Lysates can be aliquoted and snap frozen in liquid nitrogen before storing at -80°C. Avoid freeze/thaw.

### Inosine Standard:

1. Dilute Inosine Standard to 1 mM by adding 10 µl of Inosine Standard/10 mM Inosine Standard to 90 µl ADA Assay Buffer.
2. Further dilute the Inosine Standard to 10 µM by adding 10 µl of 1 mM Inosine to 990 µl ADA Assay Buffer.
3. Add 0, 2, 4, 6, 8 and 10 µl of diluted 10 µM Inosine Standard into a series of wells in a 96-well plate to generate 0, 20, 40, 60, 80 and 100 pmol/well Inosine Standard.
4. Adjust the volume to 50 µl/well with ADA Assay Buffer.

### Adenosine Deaminase Activity Assay:

1. Add 2-50 µl of sample into desired well(s) in 96-well plate. For Positive Control, dilute the Positive Control 1:10 into ADA Assay Buffer and add 1-2 µl into desired well(s).
2. Adjust the volume of sample and Positive Control to 50 µl/well with ADA Assay Buffer.
3. Add 50 µl ADA Assay Buffer to one well as reagent Background Control.

### Δ Notes:

- a. For unknown samples, we suggest doing pilot experiment and testing several doses to ensure the readings are within the Standard Curve range.
- b. Small molecules such as adenosine, inosine, xanthine, and hypoxanthine in the samples will contribute to the background. Remove these molecules by passing through a desalting column or by buffer exchange using a 10 kDa spin column. Use this modified sample for the assay. **Optional:** Prepare a parallel sample well as sample background control to ensure that these small molecules are removed by either using a desalting column or spin column.

### Reaction Mix:

1. Prepare enough reagents for the number of assays to be performed. Make 50  $\mu\text{L}$  of Reaction Mix and Background Control Mix containing:

Item	Reaction Mix	Background Control Mix
ADA Assay Buffer	40 $\mu\text{L}$	45 $\mu\text{L}$
ADA Converter	2 $\mu\text{L}$	2 $\mu\text{L}$
Converter Enzyme VIII/ADA Developer	2 $\mu\text{L}$	2 $\mu\text{L}$
OxiRed Probe/ADA Probe	1 $\mu\text{L}$	1 $\mu\text{L}$
ADA Substrate	5 $\mu\text{L}$	-

2. Add 50  $\mu\text{L}$  of Reaction Mix into each sample, reagent background control and Positive Control wells and 50  $\mu\text{L}$  of Background Control mix to Standards and sample background control well(s). Mix well.

### Measurement

Measure fluorescence (Ex/Em = 535/587 nm) in kinetic mode for at least 30 min. at 37°C. Choose two time points ( $T_1$  &  $T_2$ ) in linear range (can be as short as 2 min) of plot and obtain corresponding RFU for sample ( $\text{RFU}_{S1}$  and  $\text{RFU}_{S2}$ ) and reagent background control ( $\text{RFU}_{BG1}$  and  $\text{RFU}_{BG2}$ ). Read the Inosine Standard Curve along with the samples.

### Calculations

1. Subtract 0 Standard reading from all Standard Readings. Plot the Inosine Standard Curve.
2. Subtract reagent background control reading from sample reading.
3. Apply the  $\Delta\text{RFU}$  [ $(\text{RFU}_{S2} - \text{RFU}_{BG2}) - (\text{RFU}_{S1} - \text{RFU}_{BG1})$ ] to the Standard Curve to get B pmol of Inosine generated by the sample during the reaction time ( $\Delta T = T_2 - T_1$ ).

**Δ Note:** Sample background control reading should be less than reagent background control reading. We recommend removing the small molecules again using desalting column or a 10 kDa spin column if sample background control reading is higher than reagent background control.

$$\text{Sample ADA Activity} = \frac{B}{\Delta T \times \mu\text{g of protein}} \times \text{DF} = \text{pmol/min}/\mu\text{g} = \mu\text{U}/\mu\text{g}$$

Where: **B** = Inosine amount from the Standard Curve (pmol)  
 **$\Delta T$**  = Reaction time (min)  
 **$\mu\text{g of protein}$**  = Amount of protein/well ( $\mu\text{g}$ )  
**DF** = Dilution factor of the sample

### 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)