

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

For the detection of ADA activity.  
For research use only - not intended for diagnostic use.

Adenosine Deaminase (ADA) Activity Assay Kit (Fluorometric) (ab204695) is an assay where inosine formed from the breakdown of adenosine is detected via a multi-step reaction, resulting in the formation of an intermediate that reacts with the ADA Probe to generate a fluorescent product that can be easily quantified at Ex/Em = 535/587 nm. The kit measures total Activity of Adenosine Deaminase with limit of quantification of 5  $\mu$ U recombinant Adenosine Deaminase.

## 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	25 mL	-20°C or 4°C
Developer Solution K	1 vial	-20°C
Converter Mix F	1 vial	-20°C
ADA Positive Control	1 vial	-20°C
OxiRed™ Probe	0.2 mL	-20°C
ADA Substrate	500 $\mu$ L	-20°C
Inosine Standard	100 $\mu$ L	-20°C

PLEASE NOTE: Converter Mix F was previously labelled as Converter Enzyme VIII and ADA Developer, Developer Solution K as ADA Converter, and OxiRed™ Probe as ADA Probe. The mechanism of detection has not changed.

## 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:** Thaw completely, and ensure all particulates return to solution. Make 1x Assay Buffer by adding one part 10X ADA Assay Buffer to nine parts deionized water. Store at -20°C or 4°C. Bring to 37°C before use.

**Developer Solution K:** Reconstitute each with 220  $\mu$ l ADA Assay Buffer (1X) 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.

**Converter Mix F:** Reconstitute each with 440  $\mu$ l ADA Assay Buffer (1X) 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 Positive Control:** Reconstitute with 22  $\mu$ l ADA Assay Buffer (1X) 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. Dilute directly before use as directed in assay.

**OxiRed™ Probe:** Ready to use as supplied. Warm by placing in a 37°C bath for 1 – 5 minutes to thaw the DMSO solution before use. Keep at room temperature during the assay. Store at -20°C and protect from light and moisture. Once the probe is opened and thawed, it is stable for at least 3 additional freeze/thaw cycles but should be used within two months. After use, promptly re-tighten the cap to minimize adsorption of airborne moisture

**ADA Substrate and Inosine standard:** Ready to use when thawed. Aliquot and store at -20°C. Avoid repeated freeze/thaw.

## Assay Protocol

- Equilibrate all materials and prepared reagents to room temperature prior to use.
- **Δ Note:** Only use 1X ADA Assay Buffer diluted from 10X stock.

## 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  $\mu$ l 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  $\mu$ l 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:** Small molecules such as adenosine, inosine, xanthine, and hypoxanthine present in the samples will contribute to the background. The molecules can be removed by passing the sample through a desalting column (and using the retentate) or by buffer exchange using a 10 kDa spin column (ab93349). Use this modified processed sample for the assay.

**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  $\mu$ l of Inosine Standard to 90  $\mu$ l ADA Assay Buffer.
2. Further dilute the Inosine Standard to 10  $\mu$ M by adding 10  $\mu$ l of 1 mM Inosine to 990  $\mu$ l ADA Assay Buffer.
3. Add 0, 2, 4, 6, 8 and 10  $\mu$ l of diluted 10  $\mu$ M 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 and adjust the volume to 50  $\mu$ l/well with ADA Assay Buffer. **OR** build replicate standard curves using the table below:

Standard #	Volume of 10 $\mu$ M Standard ( $\mu$ L)*	Assay Buffer X ( $\mu$ L)	Final volume standard in well ( $\mu$ L)	End inosine Amount (pmoles/well)
1	0	125	50	0
2	5	120	50	20
3	10	115	50	40
4	15	110	50	60
5	20	105	50	80
6	25	100	50	100

Each dilution has enough volume to set up duplicate readings (2 x 50  $\mu$ L).

### Adenosine Deaminase Activity Assay:

1. Add 2-50  $\mu\text{L}$  of sample into desired well(s) in 96-well plate.
2. For Positive Control, Dilute stock by a factor of 100 (eg 2  $\mu\text{L}$  of ADA Positive control with 198  $\mu\text{L}$  of ADA Assay Buffer). and add 2-5  $\mu\text{L}$  into desired well(s).
3. Adjust the volume of sample and Positive Control to 50  $\mu\text{L}$ /well with ADA Assay Buffer.
4. Add 50  $\mu\text{L}$  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

### 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	41 $\mu\text{L}$	46 $\mu\text{L}$
Developer Solution K	2 $\mu\text{L}$	2 $\mu\text{L}$
Converter Mix F	1 $\mu\text{L}$	1 $\mu\text{L}$
OxiRed™	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. Average the duplicate reading for each standard and sample.
2. Subtract 0 Standard reading from all Standard Readings. Plot the Inosine Standard Curve. Plot the corrected absorbance values for each standard as a function of the final amount of inosine in pmoles.
3. Subtract reagent background control reading from sample reading.
4. 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.

$$\text{ADA activity} = \left( \frac{B}{\Delta T * V} \right) * D = \text{pmol}/(\text{min} * \mu\text{L}) = \mu\text{U}/\mu\text{L}$$

Where: **B** = Inosine amount from the Standard Curve (pmol)  
**ΔT** = Reaction time (min)  
**V** = Volume of sample added to well ( $\mu\text{L}$ ).  
**D** = Dilution factor of the sample

Version 9a, Last updated January 29, 2026

### Unit Definition:

**1 Unit ADA activity = amount of Adenosine Deaminase which generates 1.0  $\mu\text{mol}$  of inosine per minute at 37°C under the assay conditions.**

**Δ Note:** Once the activity per  $\mu\text{L}$  of sample is calculated, this can easily be converted to activity / cell or  $\mu\text{g}$  of tissue lysed. It may also be expressed in terms of units/ $\mu\text{g}$  of protein by dividing by the sample protein concentration (obtained with a BCA or similar protein quantification assay).

### Technical Support

Copyright © 2026 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)