

Product datasheet

Adenosine Deaminase (ADA) Activity Assay Kit (Colorimetric) ab211093

★★★★★ [1 Abreviews](#) [2 References](#) [3 Images](#)

Overview

Product name	Adenosine Deaminase (ADA) Activity Assay Kit (Colorimetric)
Detection method	Colorimetric
Sample type	Cell Lysate, Purified protein, Tissue Lysate
Assay type	Enzyme activity (quantitative)
Sensitivity	1000 µU
Species reactivity	Reacts with: Mammals, Other species
Product overview	Adenosine Deaminase (ADA) Activity Assay Kit (Colorimetric) (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 convertor and developer to generate uric acid that can be easily quantified at OD293 nm. The kit measures total activity of Adenosine Deaminase with limit of quantification of 1 mU recombinant Adenosine Deaminase.

Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K321 Adenosine Deaminase Activity Assay Kit (Colorimetric). K321-100 is the same size as the 100 test size of ab211093.</p> <p>Adenosine Deaminase (ADA, EC 3.5.4.4) is an enzyme that catalyses the conversion of adenosine and 2'-deoxyadenosine to inosine and 2'-deoxyinosine. Adenosine Deaminase is widely distributed in various tissues and cells. There are two isoforms, ADA1 and ADA2. ADA1 is widely expressed in most cells in the body, particularly in lymphocytes and macrophages. It is present in the cytosol, nucleus and has been found associated with DPP4/CD26 in the cell membrane. ADA2 was first found in the spleen but is predominantly found in the plasma and serum. Increased serum ADA levels are found in certain infectious diseases such as tuberculosis and various liver diseases such as acute hepatitis, alcoholic hepatic fibrosis, chronic active hepatitis to name a few. Adenosine Deaminase is also a marker for T-lymphocyte proliferation.</p>
Platform	Microplate reader

Properties

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

Components	100 tests
10X ADA Assay Buffer	1 x 25ml
96-Well UV Transparent Plate	1 unit
Development Enzyme Mix VIII	1 vial
ADA Positive Control	1 vial
ADA Substrate	1 x 500µl
Converter Enzyme VI	1 vial
Inosine Standard	1 x 100µl

Function Catalyzes the hydrolytic deamination of adenosine and 2-deoxyadenosine. Plays an important role in purine metabolism and in adenosine homeostasis. Modulates signaling by extracellular adenosine, and so contributes indirectly to cellular signaling events. Acts as a positive regulator of T-cell coactivation, by binding DPP4. Its interaction with DPP4 regulates lymphocyte-epithelial cell adhesion.

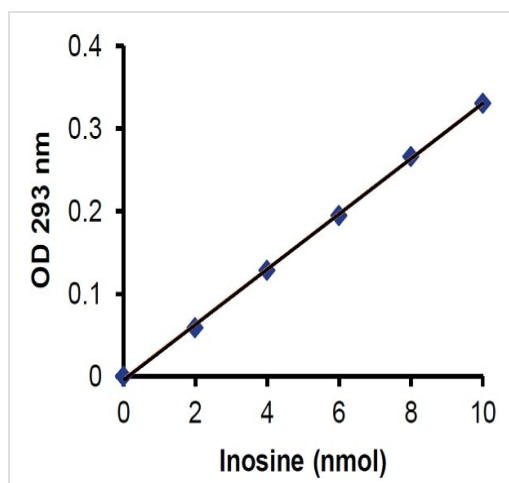
Tissue specificity Found in all tissues, occurs in large amounts in T-lymphocytes and, at the time of weaning, in gastrointestinal tissues.

Involvement in disease Defects in ADA are the cause of severe combined immunodeficiency autosomal recessive T-cell-negative/B-cell-negative/NK-cell-negative due to adenosine deaminase deficiency (ADASCID) [MIM:102700]. SCID refers to a genetically and clinically heterogeneous group of rare congenital disorders characterized by impairment of both humoral and cell-mediated immunity, leukopenia, and low or absent antibody levels. Patients with SCID present in infancy with recurrent, persistent infections by opportunistic organisms. The common characteristic of all types of SCID is absence of T-cell-mediated cellular immunity due to a defect in T-cell development. ADA-SCID is an autosomal recessive form accounting for about 50% of non-X-linked SCIDs. ADA deficiency has been diagnosed in chronically ill teenagers and adults (late or adult onset). Population and newborn screening programs have also identified several healthy individuals with normal immunity who have partial ADA deficiency.

Sequence similarities Belongs to the adenosine and AMP deaminases family.

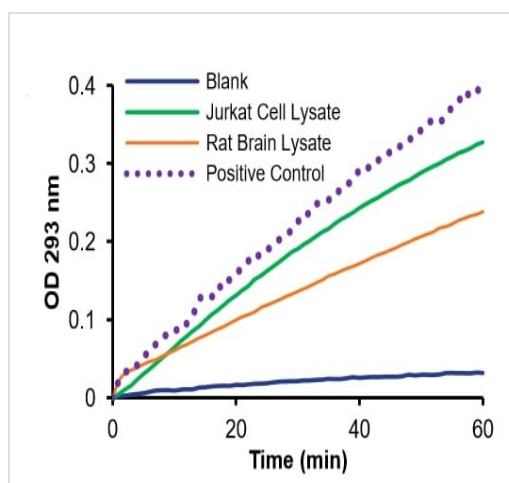
Cellular localization Cell membrane. Cell junction. Cytoplasmic vesicle lumen. Cytoplasm. Colocalized with DPP4 at the cell junction in lymphocyte-epithelial cell adhesion.

Images



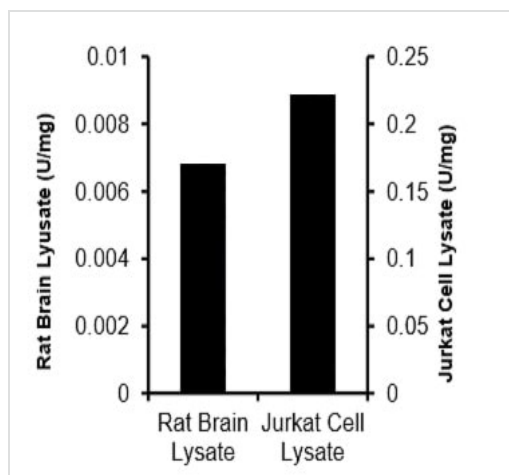
Typical inosine standard calibration curve.

Adenosine Deaminase (ADA) Activity Assay Kit
(Colorimetric) (ab211093)



Kinetic curves showing ADA activity detection in positive control (included in the kit), Jurkat cell lysate (3 μ g) and rat brain lysate (19 μ g).

Adenosine Deaminase (ADA) Activity Assay Kit
(Colorimetric) (ab211093)



ADA specific activity in rat brain lysate (19 µg) and Jurkat cell lysate (3 µg).

ADA Activity Assay Kit (Colorimetric)
(ab211093)Adenosine Deaminase (ADA) Activity
Assay Kit (Colorimetric) (ab211093)

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