

Adenosine Assay Kit (Fluorometric) ab211094

★★★★★ [3 Abreviews](#) [19 References](#) [3 Images](#)

Overview

Product name	Adenosine Assay Kit (Fluorometric)
Detection method	Fluorescent
Sample type	Urine, Plasma
Assay type	Quantitative
Assay time	0h 15m
Species reactivity	Reacts with: Mammals, Other species
Product overview	Adenosine Assay Kit (Fluorometric) (ab211094) provides a convenient method to detect adenosine in plasma and urine. In this assay, adenosine is measured using adenosine deaminase followed by a multi-step enzymatic approach resulting in the generation of an intermediate that reacts with the adenosine probe, leading to the formation of a fluorescent product. The fluorescent product can be detected at Ex/Em = 535/587 nm, and its intensity is proportional to the amount of adenosine in the sample.

The detection range of this product is 2-80 pmol of adenosine in plasma or urine.

Adenosine assay protocol summary:

- add samples and standards to wells
- add reaction mix and incubate for 15 min
- analyze with fluorescence microplate reader

Notes This product is manufactured by BioVision, an Abcam company and was previously called K327 Adenosine Assay Kit (Fluorometric). K327-100 is the same size as the 100 test size of ab211094.

Adenosine, a purine nucleoside, is present throughout the body. It plays an important role in energy transfer via the formation of ATP, ADP and AMP and in signal transduction via the formation of cAMP. Adenosine mediates its effects directly via adenosine receptors A1, A2A, A2B and A3. It regulates myocardial oxygen consumption and coronary blood flow, exerts anti-inflammatory effects throughout the body and also regulates the Renin-Angiotensin system. It also plays a role in tissue damage and repair, and cell death. Plasma adenosine levels are increased in patients with ischemic and non-ischemic heart failure.

Platform Microplate reader

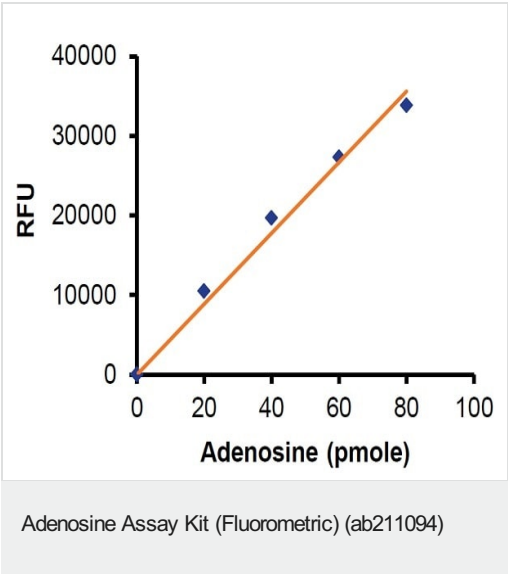
Properties

Properties

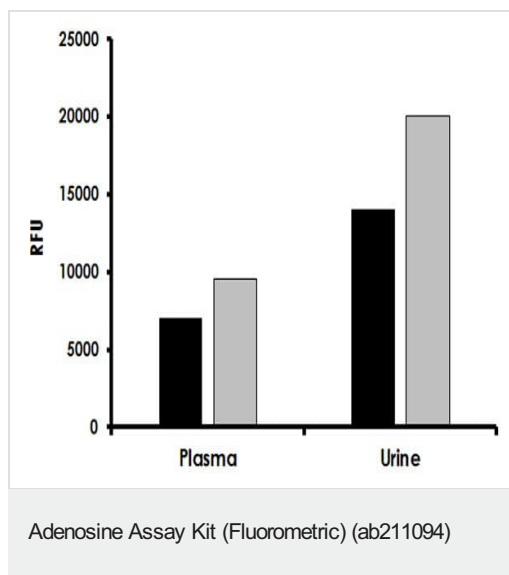
Storage instructions Store at -20°C. Please refer to protocols.

Components	100 tests
Assay Buffer X	1 x 25ml
Converter Enzyme IX	1 vial
Adenosine Deaminase I	1 vial
Adenosine Developer	1 vial
Adenosine Standard	1 x 100µl
OxiRed Probe	1 x 200µl
Xanthine Enzyme Mix	1 vial

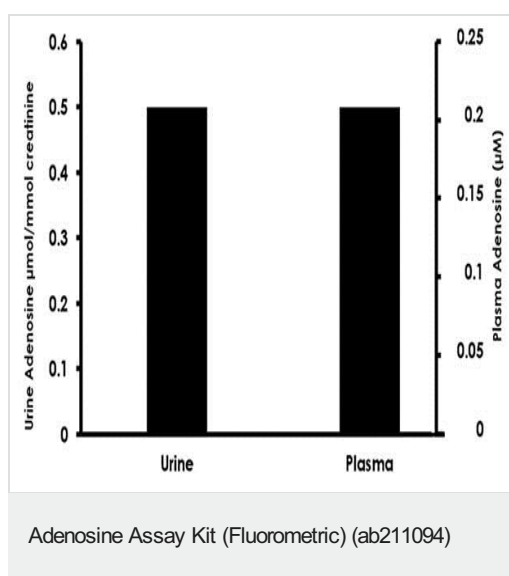
Images



Typical Adenosine standard calibration curve.



Measurement of Adenosine in pooled human plasma (20 μ L) and human urine (4 μ L of pretreated urine, 2X diluted during the pre-treatment method).



Adenosine amount in human plasma and human urine, calculated from data shown in previous figure.

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