

ADP Assay Kit (Colorimetric/Fluorometric) ab83359

27 References 4 图像

概述	
产品名称	ADP Assay试剂盒(Colorimetric/Fluorometric)
检测方法	Colorimetric/Fluorometric
样品类型	Urine, Serum, Plasma, Cell culture extracts, Other biological fluids, Tissue Extracts, Cell Lysate, Cell culture media
检测类型	Quantitative
灵敏度	> 1 μM
检测时间	1h 00m
种属反应性	与反应: Mammals, Other species
产品概述	<p>ADP Assay Kit (Colorimetric/Fluorometric) (ab83359) provides a convenient colorimetric and fluorometric method to measure ADP level.</p> <p>In the ADP assay, ADP is converted to ATP and pyruvate. The generated pyruvate can be quantified by colorimetric (ODmax = 570 nm) or fluorometric method (Ex/Em 535/587 nm).</p> <p>Conventionally, ADP levels are measured by luciferase/luciferin mediated assays after ADP is converted to ATP. However, the luciferase system is unstable and luminescence equipment is not generally available in most laboratories. In comparison, this assay is simple, sensitive, stable and high-throughput adaptable and can be used with conventional microplate readers.</p> <p>The ADP assay can detect as low as 1 μM ADP in biological samples.</p>
说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K355 ADP Colorimetric/Fluorometric Assay Kit. K355-100 is the same size as the 100 test size of ab83359.</p> <p>ADP is a product of ATP de-phosphorylation and it can be rephosphorylated to ATP. De-phosphorylation and re-phosphorylation occur via various phosphatases, phosphorylases and kinases. ADP is stored in platelets and can be released to interact with a variety of purinergic receptors. ADP levels regulate several enzymes involved in intermediary metabolism. ADP conversion to ATP primarily occurs within the mitochondrion and chloroplast although several such processes occur in the cytoplasm.</p>
平台	Microplate reader

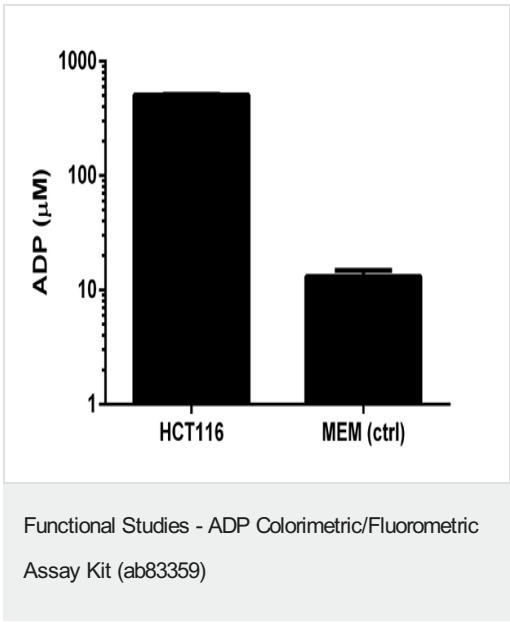
性能

存放说明 Store at -20°C. Please refer to protocols.

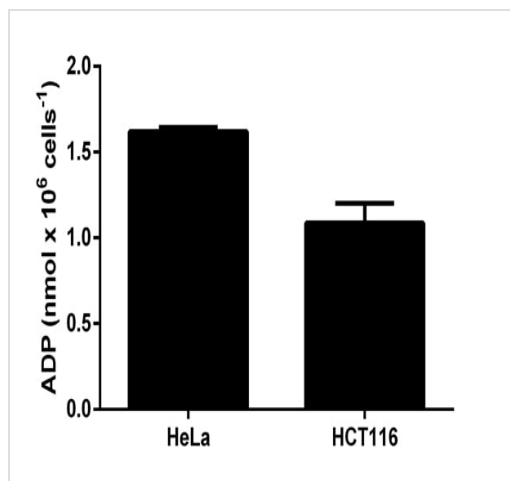
组件	100 tests
ADP Standard	1 vial
Assay Buffer IV	1 x 25ml
Converter Mix II	1 vial
Development Enzyme Mix I	1 vial
OxiRed Probe	1 x 200µl

相关性 ADP (Adenosine diphosphate) is a product of ATP dephosphorylation and it can be rephosphorylated to ATP. Dephosphorylation and rephosphorylation occur via various phosphatases, phosphorylases and kinases. ADP is stored in platelets and can be released to interact with a variety of purinergic receptors. ADP levels regulate several enzymes involved in intermediary metabolism. ADP conversion to ATP primarily occurs within the mitochondrion and chloroplast although several such processes occur in the cytoplasm.

图片

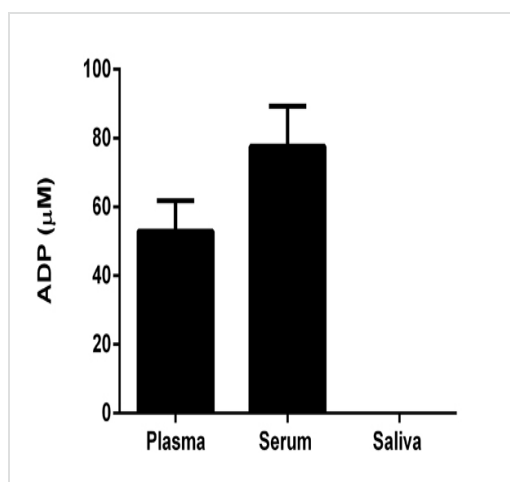


ADP levels measured fluorometrically in cell culture supernatants (background signal subtracted, mean of duplicates; +/- SD).



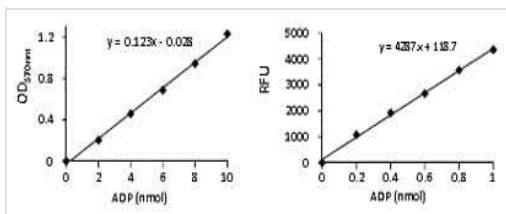
ADP levels measured fluorometrically in cell lysates (background signal subtracted, mean of duplicates; +/- SD).

Functional Studies - ADP Colorimetric/Fluorometric Assay Kit (ab83359)



ADP levels colorimetrically measured in biological fluids (rat plasma and serum, human saliva); background signal subtracted, mean of duplicates; +/- SD.

Functional Studies - ADP Colorimetric/Fluorometric Assay Kit (ab83359)



Colorimetric and Fluorometric examples of ADP Standard curve.

Functional Studies - ADP Colorimetric/Fluorometric
Assay Kit (ab83359)

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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors