

Oxaloacetate Assay Kit ab83428

9 References 3 图像

概述

产品名称	Oxaloacetate Assay试剂盒
检测方法	Colorimetric/Fluorometric
样品类型	Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids, Tissue Extracts
检测类型	Quantitative
灵敏度	> 2 μM
范围	2 μM - 200 μM
检测时间	0h 40m
产品概述	Abcam's Oxaloacetate Assay Kit provides a simple, sensitive and rapid means of quantifying OAA in a variety of samples. In the assay, OAA is converted to pyruvate which is utilized to convert a nearly colorless probe to an intensely colored (λmax= 570nm) and fluorescent (Ex/Em = 535/587nm) product. The Oxaloacetate Assay Kit can detect 0.1-10nmol (2-200 μM) of OAA.

	Visit our <a href="#">FAQs page</a> for tips and troubleshooting.
说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K659 Oxaloacetate Colorimetric/Fluorometric Assay Kit. K659-100 is the same size as the 100 test size of ab83428.</p> <p>Oxaloacetate (OAA, HOOC-CO-CH<sub>2</sub>-COOH) is a TCA cycle intermediate. It precedes citrate which is formed by the transfer of an acetyl group to OAA. OAA is formed by the deamidation of aspartate or condensation of CO<sub>2</sub> with pyruvate or PEP. Since mammals do not possess the enzymatic machinery to form TCA cycle intermediates from acetyl CoA, OAA is one of the anaplerotic entry points via pyruvate and pyruvate carboxykinase</p>

平台	Microplate reader
----	-------------------

性能

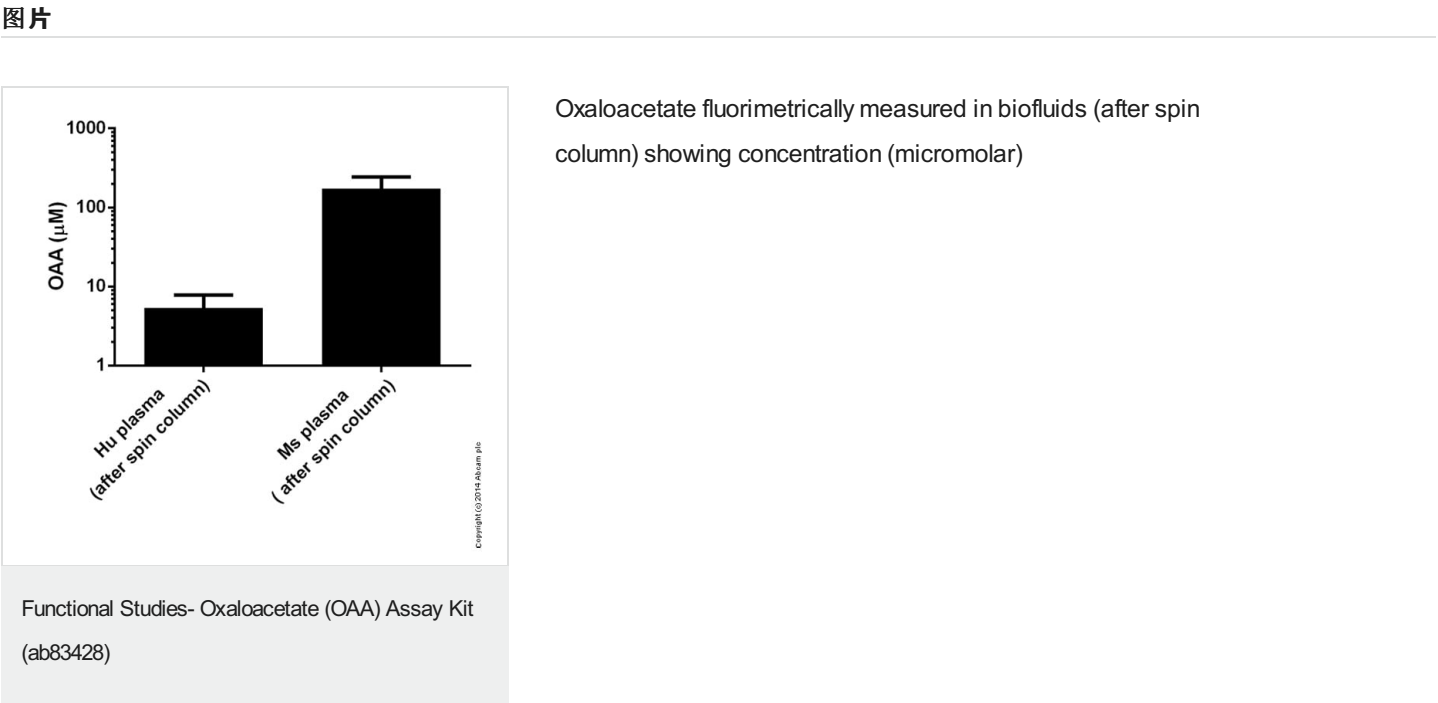
存放说明	Store at -20°C. Please refer to protocols.
------	--

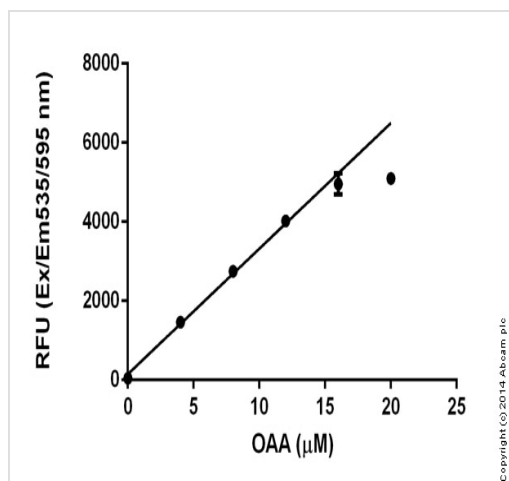
组件	100 tests
Developer	1 vial

组件	100 tests
OAA Assay Buffer	1 x 25ml
OAA Enzyme Mix	1 vial
OAA Probe (in DMSO)	1 x 200μl
OAA Standard (10 μmol)	1 vial

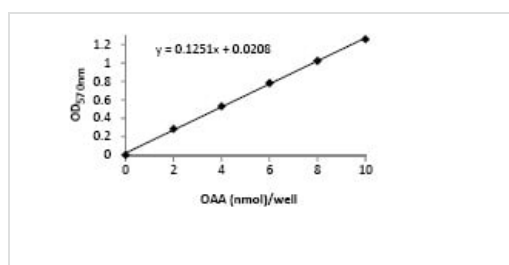
相关性

Oxaloacetate (OAA,  $\text{HOOC-CO-CH}_2\text{-COOH}$ ) is a TCA cycle intermediate. It precedes citrate which is formed by the transfer of an acetyl group to OAA. OAA is formed by the deamidation of aspartate or condensation of  $\text{CO}_2$  with pyruvate or PEP. Since mammals do not possess the enzymatic machinery to form TCA cycle intermediates from acetyl CoA, OAA is one of the anaplerotic entry points via pyruvate and pyruvate carboxykinase.





Functional Studies- Oxaloacetate (OAA) Assay Kit  
(ab83428)



Functional Studies - Oxaloacetate (OAA) Assay Kit  
(ab83428)

**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