

Product datasheet

Ascorbic Acid Assay Kit ab65346

2 图像

概述

产品名称	Ascorbic Acid Assay试剂盒
样品类型	Cell culture supernatant, Urine, Other biological fluids, Tissue Extracts
检测类型	Quantitative
范围	0.01 nmol/well - 10 nmol/well
检测时间	0h 30m
产品概述	<p>Abcam's Ascorbic Acid Assay Kit provides a rapid, simple, and sensitive means of detecting ascorbic acid in various biological samples. In this assay, our proprietary catalyst oxidizes ascorbic acid to produce a product that interacts with the ascorbic acid probe, generating color and fluorescence. Ascorbic acid can be easily determined by either colorimetric (spectrophotometry at $\lambda = 570$ nm) or fluorometric (Ex/Em = 535/587 nm) methods. The assay can detect 0.01-10 nmol of ascorbic acid per assay in various samples.</p> <p>Visit our FAQs page for tips and troubleshooting.</p>

说明	<p>Ascorbic Acid (Vitamin C) plays an important role in many biological processes. It is a potent anti-oxidant, anti-inflammatory, anti-viral agent, and an immune stimulant and is present in a wide variety of foods and biological specimens. It is important to be able to monitor ascorbic acid content in these different samples.</p>
----	--

性能

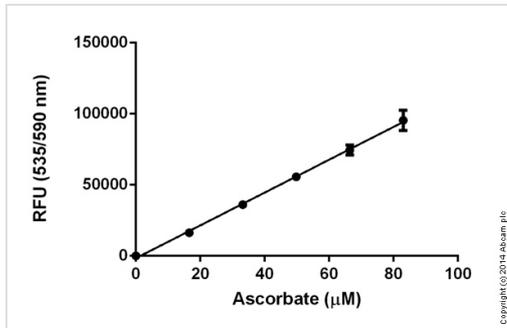
存放说明 Store at -20°C. Please refer to protocols.

组件	标识符	100 tests
Ascorbic Acid Assay Buffer	WM	1 x 25ml
Ascorbic Acid Enzyme Mix (lyophilized)	Green	1 vial
Ascorbic Acid Probe (DMSO)	Red	1 x 0.2ml
Ascorbic Acid Standard (20 μ mole)	Yellow	1 vial
Catalyst	Blue	1 x 500 μ l

相关性	<p>Ascorbic Acid (C₆H₈O₆) is an essential nutrient in human diets, and is necessary to maintain connective tissue and bone. The biologically active form of ascorbic acid is vitamic C, a water</p>
-----	--

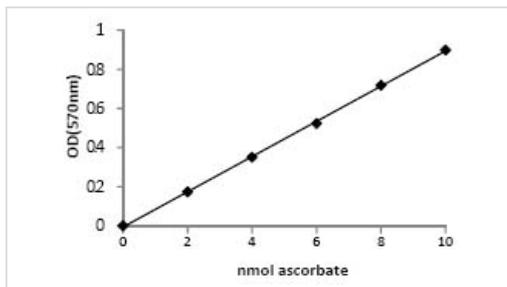
soluble vitamin which is the L-enantiomer form of ascorbic acid. Humans have lost the ability to synthesize ascorbic acid and must obtain in the diet. Ascorbic acid is an electron donor for enzymes involved in collagen hydroxylation, biosynthesis of carnitine and norepinephrine, tyrosine metabolism and amidation of peptide hormones. The ability of vitamin C to donate electrons also makes it a potent water-soluble anti-oxidant that readily scavenges free radicals.

图片



Standard curve (fluorimetric) : mean of duplicates (+/-SD) with background readings subtracted

Functional studies - Ascorbic Acid Assay Kit
(ab65346)



Ascorbic Acid Standard Curve.

Functional Studies - Ascorbic Acid Assay Kit
(ab65346)

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

Our Promise 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