

# Total Antioxidant Capacity Assay Kit ab65329

★★★★★ [3 Abreviews](#) [113 References](#) [5 图像](#)

### 概述

产品名称	Total Antioxidant Capacity Assay试剂盒
检测方法	Colorimetric
样品类型	Urine, Serum, Plasma, Other biological fluids, Tissue Extracts, Cell Lysate, Cell culture media
检测类型	Quantitative
检测时间	2h 00m
产品概述	Total Antioxidant Capacity Assay Kit ab65329 can measure either the combination of both small molecule antioxidants and proteins, or small molecules alone in the presence of our proprietary Protein Mask.

In the total antioxidant capacity assay protocol, the  $\text{Cu}^{2+}$  ion is converted to  $\text{Cu}^+$  by both small molecule and protein antioxidants. The Protein Mask prevents  $\text{Cu}^{2+}$  reduction by proteins, enabling the analysis of only the small molecule antioxidants. The reduced  $\text{Cu}^+$  ion is chelated with a colorimetric probe giving a broad absorbance peak around 570 nm, proportional to the total antioxidant capacity.

Total antioxidant capacity assay protocol summary:

- add protein mask to samples if only measuring small molecule total antioxidant capacity
- add samples and standards to wells
- add  $\text{Cu}^{2+}$  solution and incubate for 90 min at room temp
- analyze with microplate reader

### 说明

This product is manufactured by BioVision, an Abcam company and was previously called K274 Total Antioxidant Capacity (TAC) Colorimetric Assay Kit. K274-100 is the same size as the 100 test size of ab65329.

Antioxidants play an important role in preventing the formation of, and scavenging of, free radicals and other oxidizing species. There are three categories of antioxidant species: enzyme systems (GSH reductase, catalase, peroxidase, etc.), small molecules (ascorbate, uric acid, GSH, vitamin E, etc.) and proteins (albumin, transferrin, etc.).

Trolox is used to standardize antioxidants, with all other antioxidants being measured in Trolox equivalents. Measurement of the combined non-enzymatic antioxidant capacity of biological fluids and other samples provides an indication of the overall capability to counteract reactive oxygen species (ROS), resist oxidative damage and combat oxidative stress-related diseases.

### Related products

Review the [oxidative stress marker and assay guide](#) to learn about more assays for oxidative stress.

## 平台

Microplate reader

## 性能

## 存放说明

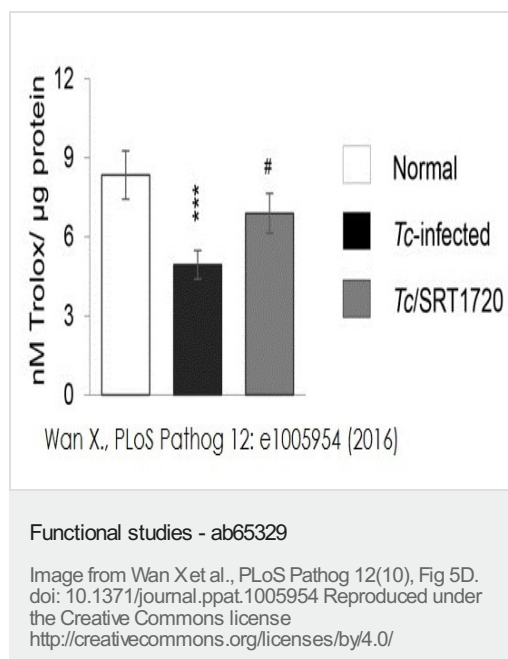
Store at +4°C. Please refer to protocols.

组件	100 tests
Assay Buffer XXIV	1 x 10ml
Copper Ion (2+) Reagent	1 x 0.2ml
Protein Mask	1 x 10ml
Trolox Standard	1 vial

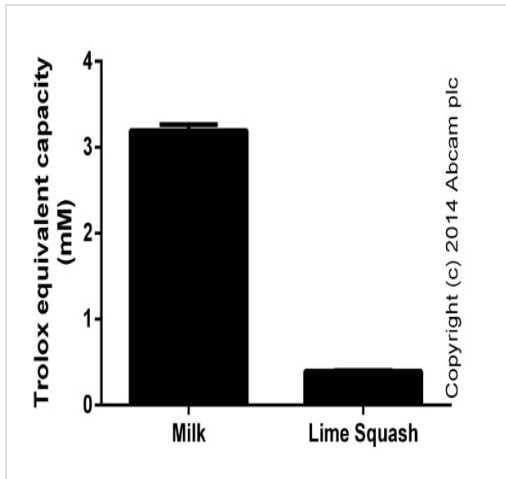
## 相关性

Antioxidants play an important role in preventing the formation of and scavenging of free radicals and other potentially toxic oxidizing species. There are three categories of antioxidant species: enzyme systems (GSH reductase, catalase, peroxidase, etc.), small molecules (ascorbate, uric acid, GSH, vitamin E, etc.) and proteins (albumin, transferrin, etc.). Different antioxidants vary in their reducing power, and cooperation of all different antioxidants provides greater protection against reactive oxygen or nitrogen radicals than any single compound alone. Therefore, the overall total antioxidant capacity may give more relevant biological information compared to that obtained by the measurement of individual components, as it considers the cumulative effect of all antioxidants present.

## 图片

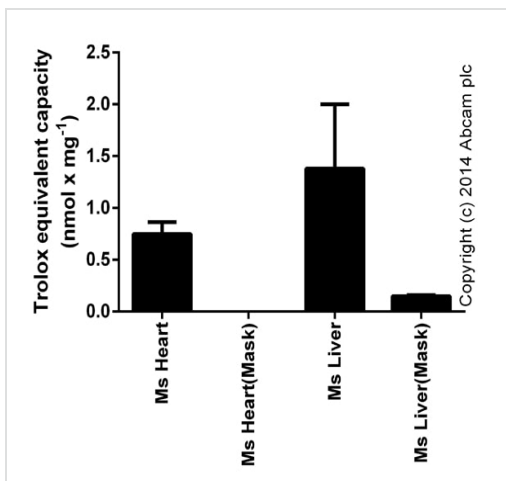


Total antioxidant capacity was determined in mouse cardiac homogenates using Total antioxidant capacity assay kit (ab65329).



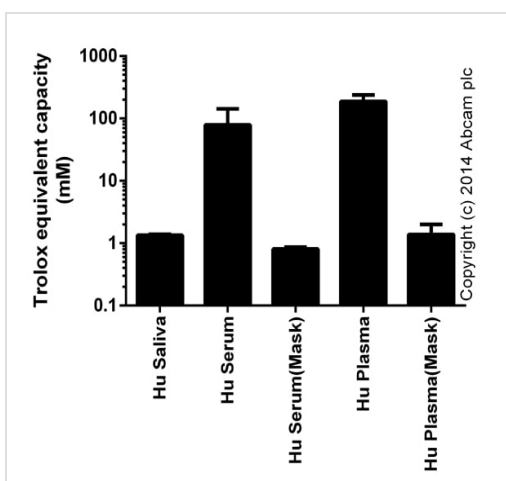
Trolox equivalent capacity measured in milk and concentrated squash. Background signal subtracted, duplicates; +/- SD

Functional assays: TAC (ab65329)



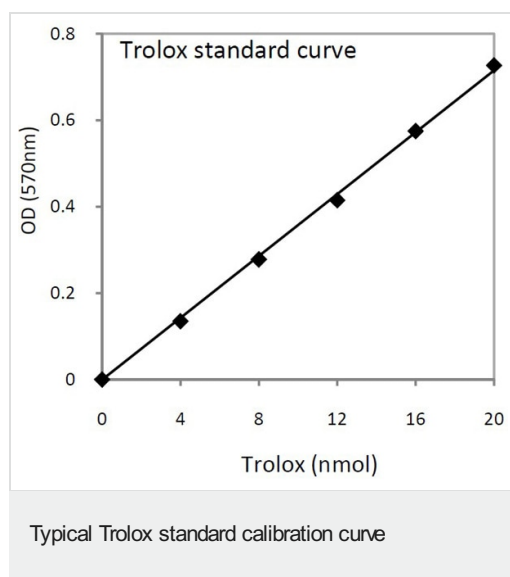
Trolox equivalent capacity measured in mouse tissue lysates, showing quantity (nmol) per mg of extracted protein. Results following blocking of protein activity is shown (Mask). (Duplicates; +/- SD).

Functional assays: TAC (ab65329)



Trolox equivalent capacity measured in biological fluids. Results following blocking of protein activity is shown (Mask). Background signal subtracted, duplicates; +/- SD.

Functional assays: TAC (ab65329)



**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