

Catalase Activity Assay Kit (Colorimetric/Fluorometric) ab83464

★★★★★ [3 Abreviews](#) [84 References](#) [2 图像](#)

概述

产品名称	Catalase Activity Assay试剂盒(Colorimetric/Fluorometric)
检测方法	Colorimetric/Fluorometric
样品类型	Urine, Serum, Plasma, Cell Lysate, Tissue Lysate
检测类型	Enzyme activity (quantitative)
检测时间	0h 40m
种属反应性	与反应: Mammals
产品概述	<p>Catalase Activity Assay Kit (Colorimetric/Fluorometric) (ab83464) is a highly sensitive, simple and direct assay for measuring catalase activity in a variety of biological samples such as cell and tissue lysates or biological fluids.</p> <p>In the catalase activity assay protocol, the catalase present in the sample reacts with hydrogen peroxide (H₂O₂) to produce water and oxygen. The unconverted H₂O₂ reacts with probe to produce a product that can be measured colorimetrically at OD 570 nm or fluorometrically at Ex/Em = 535/587 nm. Therefore, the catalase activity present in the sample is reversely proportional to the signal obtained. The kit can detect as little as 1 μU of catalase activity.</p> <p>Catalase activity assay protocol summary:</p> <ul style="list-style-type: none"> - add samples and standards to wells - add stop solution into sample control wells and incubate at 25°C for 5 min - add H₂O₂ solution into wells and incubate for 30 min at 25°C - add stop solution - add developer mix and incubate for 10 min at 25°C - analyze with microplate reader
说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K773 Catalase Activity Colorimetric/Fluorometric Assay Kit. K773-100 is the same size as the 100 test size of ab83464.</p>
平台	Microplate reader

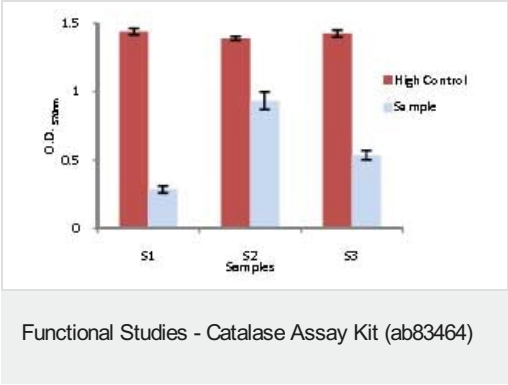
性能

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

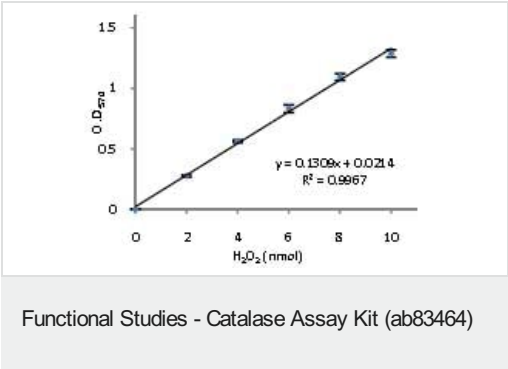
组件	100 tests	2000 tests	100 tests
Assay Buffer VIII	1 x 25ml	20 x 25ml	1 x 25ml
Catalase Positive Control	1 x 20µl	20 x 20µl	1 x 2µl
Developer Solution V	1 vial	20 vials	1 vial
Hydrogen Peroxide Solution II	1 x 25µl	20 x 25µl	1 x 25µl
OxiRed Probe	1 x 200µl	20 x 200µl	1 x 200µl
Stop Solution III	1 x 1ml	20 x 1ml	1 x 1ml

功能	Occurs in almost all aerobically respiring organisms and serves to protect cells from the toxic effects of hydrogen peroxide. Promotes growth of cells including T-cells, B-cells, myeloid leukemia cells, melanoma cells, mastocytoma cells and normal and transformed fibroblast cells.
疾病相关	Defects in CAT are the cause of acatalasia (ACATLAS) [MIM:115500]; also known as acatalasemia. This disease is characterized by absence of catalase activity in red cells and is often associated with ulcerating oral lesions.
序列相似性	Belongs to the catalase family.
翻译后修饰	The N-terminus is blocked.
细胞定位	Peroxisome.

图片



Sample tests using ab83464.



H₂O₂ standard curve using ab83464.

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