

# Complex IV Human Specific Activity Microplate Assay Kit ab109910

**11 References**   **3 图像**

### 概述

产品名称	Complex IV人Specific Activity Microplate Assay试剂盒
检测方法	Colorimetric
样品类型	Cell culture extracts, Tissue
检测类型	Enzyme activity
种属反应性	与反应: Human
产品概述	This rapid multiplexing microplate kit is used to determine both the activity and quantity of cytochrome c oxidase in a human sample with greater speed and simplicity. The ratio of the two is a measure of specific activity. The COX enzyme is immunocaptured within the wells of the microplate and the activity is determined by following the oxidation of reduced Cytochrome c colorimetrically by the absorbance change at 550 nm. Subsequently, in these same wells the quantity of the enzyme can then be determined by adding a COX specific detector antibody and the appropriate alkaline phosphatase conjugated antibody; this phosphatase changes a substrate from colorless to yellow at 405 nm. This reaction takes place in a time dependent manner proportional to the amount of protein captured in the wells. Included in this kit for performance of the multiplexing assay are buffers, detergent, substrates, and 96-well microplate with monoclonal antibody pre-bound to the wells of the plate, allowing for a stream-lined assay.

### 说明

#### Range of complex IV / cytochrome c oxidase assay kits

Biochemical assay - [ab239711](#)

Immunocapture with biochemical assay (plate-based)\*\*\* - [ab109911 \(rodent\)](#) and [ab109909 \(human\)](#)

\*\*\* Most popular assay format

Immunocapture with biochemical assay (dipstick) - [ab109878 \(rodent\)](#) and [ab109876 \(human\)](#)

Immunocapture with biochemical assay and ELISA - [ab109910 \(human\)](#) (this kit)

ELISA - [ab179880 \(human\)](#)

#### Other related products

Review the [mitochondrial assay guide](#), or the full [metabolism assay guide](#) to learn about more assays for metabolites, metabolic enzymes, mitochondrial function, and oxidative stress, and also how to assay metabolic function in live cells using your plate reader.

### 平台

Microplate reader

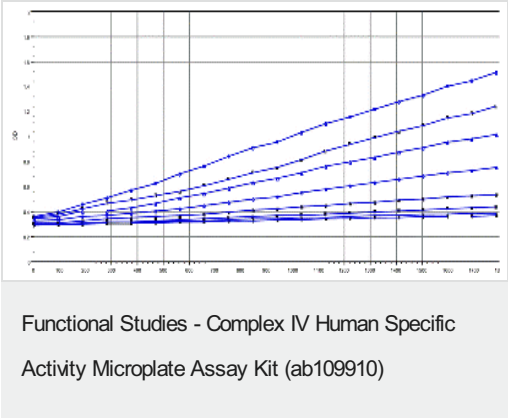
性能

存放说明

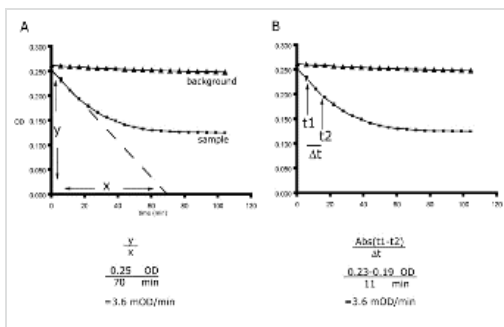
Please refer to protocols.

组件	标识符	96 tests
20X Buffer		1 x 15ml
96-well microplate (12 strips)		1 unit
AP Development Reagent		1 x 400µl
AP label 2500X	Tube B	1 x 12µl
Detector antibody	Tube A	1 x 1ml
Detergent		1 x 1ml
Development buffer	Tube 3	1 x 10ml
Reagent C (Reduced Cytochrome c)		1 x 1ml
Wash Buffer	Tube 2	1 x 2ml

图片

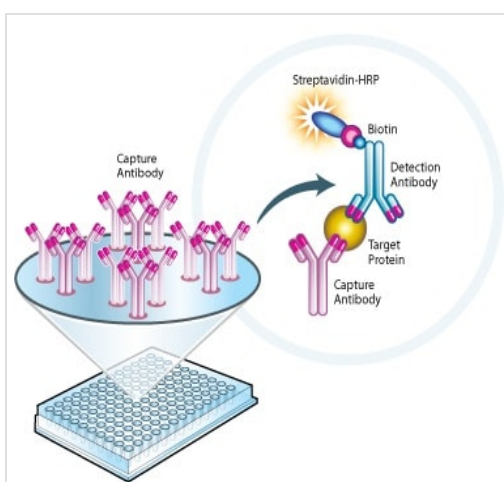


Functional study using Complex IV Human Specific Activity Microplate Assay Kit (ab109910).



Functional study using Complex IV Human Specific Activity Microplate Assay Kit (ab109910).

Functional Studies - Complex IV Human Specific Activity Microplate Assay Kit (ab109910)



Principle of Sandwich ELISA.

Sandwich ELISA - Complex IV Human Specific Activity Microplate Assay Kit (ab109910)

**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