

Product datasheet

Cytochrome c Profiling ELISA Kit [ab110172](#)

★★★★☆ 2 Abreviews 2 References 1 Image

Overview

Product name	Cytochrome c Profiling ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture extracts, Tissue
Assay type	Sandwich (qualitative)
Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Mouse, Rat, Cow, Human
Product overview	Abcam's Cytochrome C in vitro ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the measurement of Cytochrome C in Human, mouse, rat and bovine samples of purified mitochondria, tissue and cell extracts.

Abcam's Cytochrome C ELISA kit is used to determine the amount of the cytochrome c in a sample. Cytochrome c is immunocaptured within the wells and the amount is determined by adding a cytochrome c specific antibody conjugated with horseradish peroxidase (HRP). This peroxidase changes the substrate from colorless to blue. This reaction takes place in a time dependent manner proportional to the amount of protein captured in the wells. The rate of development of blue color can be followed at 600 nm or the reaction can be stopped, at a user defined time, by the addition of 100 µL of 1.5 N HCl to each well (not supplied) and read at 450 nm.

Notes	<p>The permeabilization of the mitochondrial outer membrane and the consequent release of cytochrome c and other apoptogenic proteins from the mitochondrial intermembrane space into cytoplasm is considered hallmark of many apoptotic pathways leading to cell death. Therefore assaying these proteins in mitochondrial and cytoplasmic fractions is a prime interest of many researchers. Most of the current biochemical methods of quantification of cytoplasmic and mitochondrial cytochrome c involve mechanical disruption of cells to obtain mitochondria-enriched and cytosolic fractions. These methods are time-consuming and limited to small number of samples.</p> <p>This cytochrome c detection assay was developed in conjunction with Abcam's ab109719 Cell Fractionation Kit - Standard and ab109718 Cell Fractionation Kit - HT, and the results were validated with ab110417 ApoTrack™ Cytochrome c Apoptosis Cocktail and with ab110415 ApoTrack™ Apoptosis Detection Kit for Immunocytochemistry.</p>
--------------	---

Tested applications	Suitable for: Sandwich ELISA
Platform	Microplate

Properties

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

Components	Identifier	96 tests
10X Blocking Buffer		1 x 15ml
20X Buffer	Tube 1	1 x 15ml
20X Detector antibody	Tube A	1 x 1ml
20X HRP label	Tube B	1 x 1ml
20X Wash Buffer	Tube 3	1 x 15ml
96-well pre-coated microplate		1 unit
Detergent		1 x 1ml
HRP/TMB Development Solution		1 x 20ml

Relevance Cytochrome c plays an important role in apoptosis. The protein is located in the space between the inner and outer mitochondrial membranes. An apoptotic stimulus triggers the release of cytochrome c from the mitochondria into cytosol where it binds to Apaf-1. The cytochrome c/Apaf-1 complex activates caspase-9, which then activates caspase-3 and other downstream caspases.

Cellular localization Mitochondrial

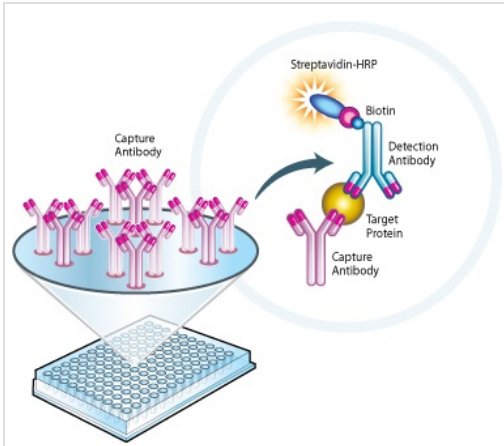
Applications

Our [Abpromise guarantee](#) covers the use of **ab110172** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
Sandwich ELISA		Use at an assay dependent dilution.

Images



Sandwich ELISA - Cytochrome c Protein Quantity
Microplate Assay Kit (ab110172)

Principle of Sandwich ELISA used in Cytochrome c Protein Quantity Microplate Assay Kit (ab110172). Substrate plus HRP or AP conjugated to streptavidin provide a colorimetric signal that is readable by any plate readers capable of standard ELISA absorbance measurements.

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

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.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors