

MitoTox™ Complex IV OXPHOS Activity Assay Kit ab109906

[6 References](#) [1 图像](#)

概述

产品名称	MitoTox™ Complex IV OXPHOS Activity Assay试剂盒
检测方法	Colorimetric
检测类型	Direct
检测时间	5h 00m
种属反应性	与反应: Cow, Human
产品概述	<p>MitoTox™ Complex IV OXPHOS Activity Assay Kit (ab109906) is designed for testing the direct inhibitory effect of compounds on Complex IV activity in only 5 hours. Complex IV extracted from the provided bovine heart mitochondria (a rich source of Complex IV) is immunocaptured by specific antibodies on the plate. Complex IV activity can be observed as decrease in absorbance at OD 550 nm. The intra-assay and inter-assay variation of this assay are both <10%.</p> <p>Inhibitory effects of compounds on Complex IV activity can be tested in two different ways: 1. Screening format, where up to 23 compounds can be tested at a single concentration in triplicate; 2. Dose response (IC₅₀) format, where two compounds known to affect Complex IV activity can be tested at 11 different data points in triplicate.</p> <p>Testing for mitochondrial function has become a key aspect of drug discovery. Mitochondria can be affected by drug treatment, resulting into cardio- and hepatotoxic side effects that can lead to drug withdrawal from the market. Therefore, there is increasing emphasis on testing the impact on mitochondria early on in the drug development process to reduce failure rates during preclinical and clinical phases.</p>
说明	<p>Store Bovine Heart Mitochondria and Reagent C at -80°C. Store all other components at 4°C.</p> <p>Related products</p> <p>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.</p>
平台	Microplate reader

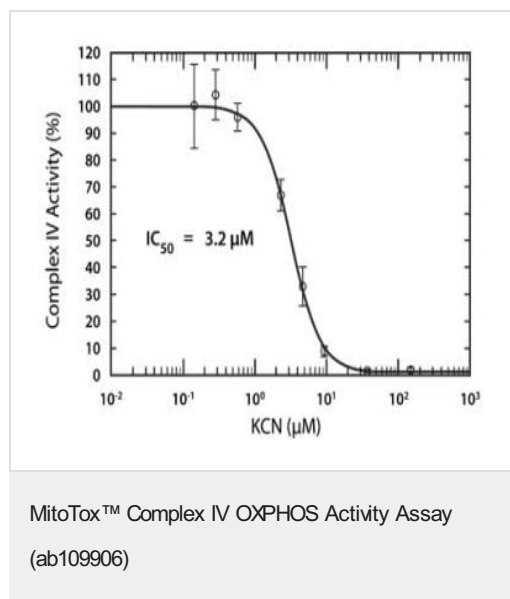
性能

存放说明

Please refer to protocols.

组件	96 tests
12-channel reagent reservoirs	2 units
1X Blocking Buffer	1 x 70ml
1X Complex IV Mito Buffer	1 x 110ml
Bovine Heart Mitochondria	1 x 90 μ l
Detergent	1 x 20 μ l
Blocking Powder	1 x 1.8g
Pre-coated 96-well microplate	1 unit
Reagent C (Reduced Cytochrome c)	2 x 1ml

图片



Typical dose response curve for KCN. Assay was performed following the Dose Response Assay Procedure using KCN, a well known Complex IV inhibitor. KCN was prepared in 0.1 M NaOH to generate a 100 mM stock. Starting with a 100 μ M final concentration in well, 1:10 serial dilutions of KCN were generated.

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