

# MitoTox™ Complex I OXPHOS Activity Assay Kit ab109903

**13 References** [画像数 1](#)

### 製品の概要

製品名	MitoTox™ Complex I OXPHOS Activity Assay Kit
検出方法	Colorimetric
サンプルの種類	Inhibitor compounds
アッセイタイプ	Quantitative
全工程の試験時間	5h 00m
製品の概要	<p>MitoTox™ Complex I OXPHOS Activity Assay (ab109903) is designed for testing the direct inhibitory effect of compounds on Complex I activity in only 5 hours. Complex I extracted from the provided bovine heart mitochondria (a rich source of Complex I) is immunocaptured by specific antibodies on the plate. Complex I activity can be observed as decrease in absorbance at OD 340 nm. The intra-assay and inter-assay variation of this assay are both &lt; 10%.</p> <p>Inhibitory effects of compounds on Complex I 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 (IC<sub>50</sub>) format, where two compounds known to affect Complex I 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>
特記事項	<p>Store Phospholipids, Bovine heart mitochondria, Ubiquinone and Activity Buffer at -80°C. Store all other components at 4°C.</p> <p><b>Related products</b></p> <p>Review the <a href="#">mitochondrial assay guide</a>, or the full <a href="#">metabolism assay guide</a> 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>
試験プラットフォーム	Microplate reader

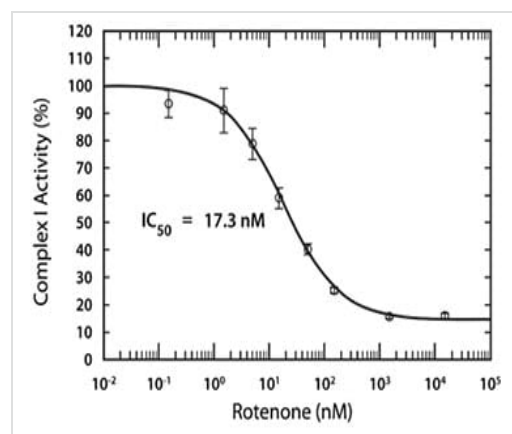
### 製品の特性

## 保存方法

Please refer to protocols.

内容	96 tests
12-channel reagent reservoirs	2 units
1X Mito Buffer	1 x 5ml
20X Wash Buffer	1 x 5ml
Bovine heart mitochondria	1 x 360µl
Complex I Activity Buffer	1 x 24ml
Detergent	1 x 100µl
Phospholipids	1 x 6ml
Pre-coated 96-well microplate	1 unit
Ubiquinone 1	1 x 60µl

## 画像



MitoTox™ Complex I OXPHOS Activity Assay  
(ab109903)

Typical dose response curve for rotenone. Assay was performed following the Dose Response Assay Procedure using rotenone, a well known Complex I inhibitor. Rotenone was prepared in DMSO to generate a 10 mM stock. Starting with a 50 µM final concentration in well, 1:10 serial dilutions of rotenone 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.co.jp/abpromise> or contact our technical team.

#### **Terms and conditions**

---

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