

A convenient, easy-to-follow shortened protocol is provided with this assay.
For a detailed protocol go to www.caymanchem.com/pdfs/700930.pdf

MitoCheck® Complex I Activity Assay Kit Short Protocol

Item No. 700930

REAGENT PREPARATION

- Mitochondrial Complex I Activity Assay Buffer - (Item No. 700931)** - Ready to use as supplied; warm to room temperature and vortex prior to use.
- NADH Assay Reagent - (Item No. 700932)** - Dissolve in 155 μ l of UltraPure water; stable on ice for three hours; store at -20°C for up to 2 weeks.
- Mitochondrial Inhibitors - (Not Supplied)**
KCN-Use extreme care when preparing the KCN reagent. Weigh 6.5 mg KCN in a ventilated hood; dissolve in 1 ml of 0.1 M NaOH to make a 500 mM stock solution; store on ice; make fresh 3 hrs before use.
Rotenone-To ensure inhibition of complex I, use concentrations ≥ 1 μ M; can be made up in DMSO or ethanol; if making up in DMSO, avoid freeze/thaws. Use appropriate PPE.
Antimycin A-To ensure inhibition of complex III, use concentrations ≥ 10 μ M; can be made up in DMSO or ethanol. Use appropriate PPE.
- Reaction Stock Solutions** - Add the following reagents into 2 polystyrene tubes (sufficient for 20 reactions).

Tube A (1 mL)	Tube B (675 mL)
910 μ l of Complex I Activity Assay Buffer	625 μ l of Complex I Activity Assay Buffer
20 μ l of 100 mM KCN (1 mM)	30 μ l of NADH Assay Reagent
50 μ l of FF-BSA Assay Reagent	20 μ l of Ubiquinone Assay Reagent
20 μ l Bovine Heart Mitochondria Assay Reagent*	

*Isolated mitochondria can settle over time; mix well before use.



Short Protocol
Item 700930

Tech Support
888.526.5351
techserv@caymanchem.com

For Research Use Only
www.caymanchem.com

MitoCheck® Complex I Activity Assay Kit Short Protocol

Item No. 700930

PERFORMING THE ASSAY

1. Add 50 µl of the contents of tube A to each well.
2. Add 20 µl of test compounds, positive control, or vehicle diluted in Assay Buffer to the appropriate wells. Allow for pre-incubation if required.
3. Add 30 µl of the contents of tube B to each well to start the reaction.
4. Immediately place plate on plate reader and measure absorbance at 340 nm (30 second intervals for 15 min @ 25°C).

CALCULATIONS

1. Plot time-dependent reaction data as absorbance (y-axis) *versus* time (x-axis).
2. To determine the reaction rate, calculate the slope for the linear portion of the curve.
3. Determine % activity using the equation below.
4. To generate a concentration response curve, plot the % activity as a function of test compound concentration.

$$\text{Complex I Activity (\%)} = \left[\frac{\text{Rate of Sample wells}}{\text{Rate of Vehicle Control}} \right] \times 100$$



Short Protocol
Item 700930

Tech Support
888.526.5351
techserv@caymanchem.com

For Research Use Only
www.caymanchem.com

07/10/2015