

A convenient, easy-to-follow shortened protocol is provided with this assay.
For a detailed protocol go to www.caymanchem.com/pdfs/700940.pdf

MitoCheck® Complex II Activity Assay Kit Short Protocol

Item No. 700940

REAGENT PREPARATION

- Mitochondrial Complex II Activity Assay Buffer - (Item No. 700941)** - Ready to use as supplied; warm to room temperature and vortex prior to use.
- 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 $\geq 1 \mu\text{M}$; 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 $\geq 10 \mu\text{M}$; can be made up in DMSO or ethanol. Use appropriate PPE.
2-Thenoyltrifluoroacetone (TTFA)-To ensure inhibition of complex II, use concentrations $\geq 1 \text{ mM}$; 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 μL)
956 μL of Complex II Activity Assay Buffer	487 μL of Complex II Activity Assay Buffer
20 μL Bovine Heart Mitochondria Assay Reagent*	8 μL of Succinate Assay Reagent
2 μL of 1 mM Rotenone	20 μL of Ubiquinone Assay Reagent
20 μL of 100 mM KCN (1 mM)	160 μL of DCPIP Assay Reagent
2 μL of 10 mM Antimycin A	

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



Short Protocol
Item 700940

Tech Support
888.526.5351
techserv@caymanchem.com

For Research Use Only
www.caymanchem.com

MitoCheck® Complex II Activity Assay Kit Short Protocol

Item No. 700940

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. Incubate five minutes at 25°C, then place plate in plate reader and measure absorbance at 600 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.

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



Short Protocol
Item 700940

Tech Support
888.526.5351
techserv@caymanchem.com

For Research Use Only
www.caymanchem.com

07/10/2015