

A convenient, easy-to-follow shortened protocol is provided with this assay.  
For a detailed protocol go to [www.caymanchem.com/pdfs/701000.pdf](http://www.caymanchem.com/pdfs/701000.pdf)

## MitoCheck® Complex V Activity Assay Kit Short Protocol

### Item No. 701000

#### REAGENT PREPARATION

1. **Mitochondrial Complex V Activity Assay Buffer - (Item No. 701001)** - Ready to use as supplied; warm to room temperature and vortex prior to use.
2. **Mitochondrial Complex V Enzyme Mix - (Item No. 701002)** - Ready to use as supplied; thaw on ice prior to use.
3. **Bovine Heart Mitochondria Assay Reagent - (Item No. 700019)** - Ready to use as supplied; thaw on ice prior to use.
4. **Mitochondrial Complex V NADH Reagent - (Item No. 701003)** - Reconstitute vial with 120 µl UltraPure water prior to use.
5. **Mitochondrial Complex V NADH Reagent - (Item No. 701003)** - Reconstitute vial with 120 µl UltraPure water prior to use
6. **Mitochondrial Inhibitors - (Not Supplied)**  
Oligomycin-Use stock at 1 mg/ml; make fresh 3 hrs before use.  
Rotenone-To ensure inhibition of complex I, use concentrations  $\geq 1$  µM; can be made up in DMSO or ethanol; if making up in DMSO, avoid freeze/thaws. Use appropriate PPE.
7. **Reaction Stock Solutions** - Add the following reagents into 2 polystyrene tubes (sufficient for 20 reactions).

Tube A (1 mL)	Tube B (675 mL)
978 µl of Complex V Activity Assay Buffer	635 µl of Complex V Assay Enzyme Mix
20 µl Bovine Heart Mitochondria Assay Reagent*	20 µl of Complex V ATP Reagent
20 µl Rotenone - not supplied (from 1 mM stock)	20 µl of Complex V NADH Reagent

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



Short Protocol  
**Item 701000**

**Tech Support**  
888.526.5351  
[techserv@caymanchem.com](mailto:techserv@caymanchem.com)

For Research Use Only  
[www.caymanchem.com](http://www.caymanchem.com)

# MitoCheck® Complex V Activity Assay Kit Short Protocol

## Item No. 701000

### PERFORMING THE ASSAY

1. Add 50 µl of the contents of tube A to each well.
2. Add 20 µl of positive control or test compounds. 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 30 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.

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



Short Protocol  
Item 701000

Tech Support  
888.526.5351  
techserv@caymanchem.com

For Research Use Only  
[www.caymanchem.com](http://www.caymanchem.com)

07/10/2015