

## Product Information

### Aconitase Activity Assay Kit

Catalog Number **MAK051**  
Storage Temperature 2–8 °C

## TECHNICAL BULLETIN

### Product Description

The aconitases are a family of iron-sulfur containing enzymes that catalyze the isomerization of citrate to isocitrate. Two aconitases, encoded by different genes, have been identified in eukaryotes. Aconitase 1 (ACO1, c-aconitase) is a cytosolic enzyme, while Aconitase 2 (ACO2, m-aconitase) is a mitochondrial enzyme that functions in the tricarboxylic acid cycle. Aconitases are reversibly inactivated by oxidative stress, and aconitase activity in cells and tissues has been used as a biomarker for oxidative damage.

The Aconitase Activity Assay kit provides a simple and direct procedure for measuring Aconitase activity in a variety of samples. Aconitase activity is determined in a coupled enzyme reaction in which citrate is converted to isocitrate by aconitase. This results in a colorimetric (450 nm) product proportional to the enzymatic activity present. One unit of aconitase is the amount of enzyme that will isomerize 1.0  $\mu$ mole of citrate to isocitrate per minute at pH 7.4 at 25 °C.

### Components

The kit is sufficient for 100 assays in 96 well plates.

|                        |        |
|------------------------|--------|
| Assay Buffer           | 25 mL  |
| Catalog Number MAK051A |        |
| Substrate              | 1 vL   |
| Catalog Number MAK051B |        |
| Developer              | 1 vL   |
| Catalog Number MAK051C |        |
| Enzyme Mix             | 0.6 mL |
| Catalog Number MAK051D |        |
| Cysteine               | 1 vL   |
| Catalog Number MAK051E |        |

(NH<sub>4</sub>)Fe(SO<sub>4</sub>)<sub>2</sub> 1 vL  
Catalog Number MAK051F

Isocitrate Standard, 100 mM 100  $\mu$ L  
Catalog Number MAK051G

### Reagents and Equipment Required but Not Provided.

- 96 well flat-bottom plate – It is recommended to use clear plates for colorimetric assays.
- Spectrophotometric multiwell plate reader

### Precautions and Disclaimer

For R&D use only. Not for drug, household, or other uses. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

### Preparation Instructions

Briefly centrifuge vials before opening. Use ultrapure water for the preparation of reagents. To maintain reagent integrity, avoid repeated freeze/thaw cycles.

Assay Buffer – Allow buffer to come to room temperature before use.

Substrate – Reconstitute with 220  $\mu$ L of water. Mix well by pipetting (don't vortex), then aliquot and store, protected from light, at 2–8 °C. Use within 2 months of reconstitution.

Developer – Reconstitute with 1.1 mL of Assay Buffer. Mix well by pipetting (don't vortex), then aliquot and store, protected from light, at 2–8 °C. Use within 2 months of reconstitution.

Aconitase Activation Solution – Reconstitute cysteine and  $(\text{NH}_4)\text{Fe}(\text{SO}_4)_2$  each with 0.5 mL of Assay Buffer. Mix well by pipetting (don't vortex), then aliquot and store, protected from light, at  $-20^\circ\text{C}$ .

To prepare Aconitase Activation Solution, mix 0.1 mL each of the reconstituted cysteine and  $(\text{NH}_4)\text{Fe}(\text{SO}_4)_2$  solutions, and mix well by pipetting. Activation solution should be **prepared fresh**. Keep on ice while in use.

### Storage/Stability

The kit is shipped on wet ice and storage at  $2-8^\circ\text{C}$ , protected from light, is recommended.

### Procedure

All samples and standards should be run in duplicate.

#### Isocitrate Standards for Colorimetric Detection

Dilute 10  $\mu\text{L}$  of the 100 mM Isocitrate Standard solution with 490  $\mu\text{L}$  of the Assay Buffer to prepare a 2 mM standard solution. Add 0, 2, 4, 6, 8, and 10  $\mu\text{L}$  of the 2 mM standard solution into a 96 well plate, generating 0 (blank), 4, 8, 12, 16, and 20 nmole/well standards. Add Assay Buffer to each well to bring the volume to 50  $\mu\text{L}$ .

#### Sample Preparation

Tissue (20–40 mg) or cells ( $1 \times 10^6$ ) can be homogenized in 100  $\mu\text{L}$  of ice-cold Assay Buffer. Centrifuge the samples at  $800 \times g$  for 10 minutes at  $4^\circ\text{C}$  to remove insoluble material. The supernatant can be used for the c-aconitase assay. For the m-aconitase assay, centrifuge the supernatant at  $20,000 \times g$  for 15 minutes at  $4^\circ\text{C}$ , collect the pellet, and dissolve in 0.1 mL cold Assay Buffer. Sonicate pellet for 20 seconds. Store samples at  $-80^\circ\text{C}$  and keep on ice while in use.

To activate samples, add 10  $\mu\text{L}$  of Aconitase Activation Solution to 100  $\mu\text{L}$  of sample and incubate on ice for 1 hour.

**Note:** For unknown samples, it is suggested to test several sample dilutions to ensure the readings are within the linear range of the standard curve.

Add 2–50  $\mu\text{L}$  of activated samples to wells. Bring samples to a final volume of 50  $\mu\text{L}$  with Assay Buffer. To control for background activity, a sample blank may be set up for each sample by omitting the Substrate (Table 1). The blank readings can be subtracted from the sample readings.

#### Assay Reaction

1. Set up the Reaction Mixes according to the scheme in Table 1. 50  $\mu\text{L}$  of the appropriate Reaction Mix is required for each reaction (well). The enzyme mix should be kept on ice while in use.

**Table 1.**  
Reaction Mixes

| Reagent      | Samples and Standards | Sample Blank     |
|--------------|-----------------------|------------------|
| Assay Buffer | 46 $\mu\text{L}$      | 48 $\mu\text{L}$ |
| Enzyme Mix   | 2 $\mu\text{L}$       | 2 $\mu\text{L}$  |
| Substrate    | 2 $\mu\text{L}$       | –                |

2. Add 50  $\mu\text{L}$  of the appropriate Reaction Mix to each of the wells. Mix well using a horizontal shaker or by pipetting. Protect the plate from light during the incubation
3. Incubate the plate at  $25^\circ\text{C}$  for 30–60 minutes.
4. Add 10  $\mu\text{L}$  of the Developer to each well, mix, and incubate at  $25^\circ\text{C}$  for 10 minutes. Measure the absorbance at 450 nm.

## Results

### Calculations

Correct for the background by subtracting the value obtained for the 0 (blank) Isocitrate standard from all sample readings. Background values can be significant and must be subtracted from all readings. Plot the Isocitrate standard curve.

Note: A new standard curve must be set up each time the assay is run.

Subtract the blank sample value from the sample readings to obtain the corrected measurement. Using the corrected measurement, the amount of isocitrate generated by the aconitase in the sample may be determined from the standard curve.

The aconitase activity of a sample may be determined by the following equation:

$$\text{Aconitase Activity (milliunit/mL)} = \frac{B \times \text{Sample Dilution Factor}}{T \times V}$$

B = Amount (nmole) of isocitrate generated

T = time reaction incubated in minutes

V = pretreated sample volume (mL) added to well

Aconitase activity is reported as nmole/min/mL = milliunit/mL. One unit of aconitase is the amount of enzyme that will isomerize 1.0  $\mu$ mole of citrate to isocitrate per minute at pH 7.4 at 25 °C.

Example:

Isocitrate (B) = 5.84 nmole

T = 45 minutes

Sample volume (V) = 0.05 mL

Sample dilution is 1

Aconitase activity is:

$$\frac{5.84 \times 1}{45 \times 0.05} = 2.59 \text{ milliunits/mL}$$

**Troubleshooting Guide**

| <b>Problem</b>                                 | <b>Possible Cause</b>                                     | <b>Suggested Solution</b>                                                                  |
|------------------------------------------------|-----------------------------------------------------------|--------------------------------------------------------------------------------------------|
| Assay not working                              | Cold assay buffer                                         | Assay Buffer must be at room temperature                                                   |
|                                                | Omission of step in procedure                             | Refer and follow Technical Bulletin precisely                                              |
|                                                | Plate reader at incorrect wavelength                      | Check filter settings of instrument                                                        |
|                                                | Type of 96 well plate used                                | For colorimetric assays, use clear plates                                                  |
| Samples with erratic readings                  | Samples prepared in different buffer                      | Use the Assay Buffer provided                                                              |
|                                                | Cell/Tissue culture samples were incompletely homogenized | Repeat the sample homogenization, increasing the length and extent of homogenization step. |
|                                                | Samples used after multiple freeze-thaw cycles            | Aliquot and freeze samples if needed to use multiple times                                 |
|                                                | Presence of interfering substance in the sample           | If possible, dilute sample further                                                         |
|                                                | Use of old or inappropriately stored samples              | Use fresh samples and store correctly until use                                            |
| Lower/higher readings in samples and standards | Improperly thawed components                              | Thaw all components completely and mix gently before use                                   |
|                                                | Use of expired kit or improperly stored reagents          | Check the expiration date and store the components appropriately                           |
|                                                | Allowing the reagents to sit for extended times on ice    | Prepare fresh Master Reaction Mix before each use                                          |
|                                                | Incorrect incubation times or temperatures                | Refer to Technical Bulletin and verify correct incubation times and temperatures           |
|                                                | Incorrect volumes used                                    | Use calibrated pipettes and aliquot correctly                                              |
| Non-linear standard curve                      | Use of partially thawed components                        | Thaw and resuspend all components before preparing the reaction mix                        |
|                                                | Pipetting errors in preparation of standards              | Avoid pipetting small volumes                                                              |
|                                                | Pipetting errors in the Reaction Mix                      | Prepare a Master Reaction Mix whenever possible                                            |
|                                                | Air bubbles formed in well                                | Pipette gently against the wall of the plate well                                          |
|                                                | Standard stock is at incorrect concentration              | Refer to the standard dilution instructions in the Technical Bulletin                      |
|                                                | Calculation errors                                        | Recheck calculations after referring to Technical Bulletin                                 |
|                                                | Substituting reagents from older kits/lots                | Use fresh components from the same kit                                                     |
| Unanticipated results                          | Samples measured at incorrect wavelength                  | Check the equipment and filter settings                                                    |
|                                                | Samples contain interfering substances                    | If possible, dilute sample further                                                         |
|                                                | Sample readings above/below the linear range              | Concentrate or dilute samples so readings are in the linear range                          |

SS,TD,MF,LS,MAM 06/19-1