

# Isocitrate Colorimetric Assay Kit

(Catalog #K656-100; 100 assays; Store Kit at -20°C)

**I. Introduction:**

Isocitric acid (HOOC-CHOH-CH (-COOH)-CH<sub>2</sub>-COOH) is an intermediate of the Krebs TCA cycle, positioned between citrate and α-ketoglutarate. It is the branch point from which the glyoxylate shunt operates in plants and lower organisms. Isocitrate is found in substantial concentrations in many fruits and vegetables as well as in foods produced from these raw materials. In the TCA cycle, isocitrate is oxidized by isocitrate dehydrogenase (IDH) to α-ketoglutarate with the generation of NAD(P)H. Loss of NAD-IDH has been implicated as a potential causative factor in retinitis pigmentosa. BioVision's Isocitrate Colorimetric Assay Kit provides a simple, sensitive and rapid means of quantifying isocitrate in a variety of samples. In this assay, isocitrate is oxidized with the generation of NADPH, which converts a nearly colorless probe to an intensely colored species with a λ<sub>max</sub> of 450 nm. The Isocitrate Assay Kit can detect 1 to 20 nmole (~0.2-5 µg) of isocitrate.

**II. Kit Contents:**

Components	K656-100	Cap Code	Part Number
Isocitrate Assay Buffer	25 ml	WM	K656-100-1
Isocitrate Enzyme Mix	200 µl	Green	K656-100-2
Developer	lyophilized	Purple	K656-100-3
Isocitrate Standard (100 mM)	100 µl	Yellow	K656-100-4

**III. Storage and Handling:**

Store kit at -20°C, protected from light. Warm Isocitrate Assay Buffer to room temperature (RT) before use. Briefly centrifuge all small vials prior to opening.

**IV. Reagent Preparation and Storage Conditions:**

**Isocitrate Enzyme Mix:** Ready to use as supplied. Use within two months.

**Developer:** Add 220 µl dH<sub>2</sub>O. Pipette up and down several times to completely dissolve (don't vortex). Store at -20°C, protected from light. Use within 2 months.

**Isocitrate Standard:** Ready to use as supplied. Keep cold while in use. Store at -20°C.

**V. Assay Protocol:**

**1. Standard Curve Preparations:**

Dilute Isocitrate Standard to 2 nmol/µl by adding 20 µl of the Isocitrate Standard to 980 µl of dH<sub>2</sub>O, mix well. Add 0, 2, 4, 6, 8, 10 µl into a series of wells of a 96 well plate. Adjust the volume to 50 µl/well with Isocitrate Assay Buffer to generate 0, 4, 8, 12, 16, 20 nmol/well of the Isocitrate Standard.

**2. Sample Preparation:**

**Tissue or Cells:** Tissue 20 mg or cells (2 x 10<sup>6</sup>) should be rapidly homogenized with 100 µl Isocitrate Assay Buffer. Centrifuge at 15,000 g for 10 min at 4°C to remove any cell debris. Enzymes in samples may interfere with the assay. We suggest deproteinizing your sample(s) using a Perchloric Acid/KOH protocol (BioVision, Cat. # K808-200) or 10 kDa molecular weight cut off spin columns (BioVision, Cat. # 1997-25). Add 1-50 µl samples into duplicate wells of a 96-well plate. Bring volume to 50 µl using Isocitrate Assay Buffer.

**Food or Beverage samples:** Most beverages can be used directly in the assay, with appropriate dilution. In general, samples should be spin filtered through a 10kDa MWCO filter such as BioVision Cat. # 1997-25. This will remove inhibitory substances, protein and most color. Solids should be processed by homogenizing 20 mg with 500 µl distilled water, with mild heating for 30 min, then centrifuge 15,000x g, 10 min, take supernatant, spin filter and dilute appropriately for the assay. For all samples, we suggest testing several doses of your samples to ensure readings are within the Standard Curve range.

**3. Reaction Mix:** Mix enough reagents for the number of samples and Standards to be performed: For each well, prepare a total 50 µl Reaction Mix containing:

Isocitrate Assay Buffer	46 µl
Isocitrate Enzyme Mix	2 µl
Developer	2 µl

\*\* NADH and NADPH can generate significant background in some instances. If interfering levels of these are suspected of being in the sample, a background control can be performed by running a parallel sample with the Isocitrate Enzyme Mix being omitted.

Add 50 µl of Reaction Mix to each well containing the Isocitrate Standard and test and background control samples.

**4. Incubate for 30 min at 37°C, protected from light.**

**5. Measure OD at 450 nm in a microplate reader.**

**6. Calculation:** Correct background by subtracting the value of the 0 Isocitrate Standard from all readings. (Note: The background reading can be significant and must be subtracted.) Plot the Standard Curve. Then apply the corrected sample readings to the Standard Curve to get Isocitrate amount in the sample wells.

The Isocitrate concentrations in the test samples:

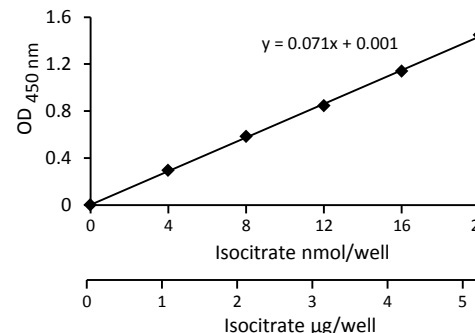
$$C = Ay/Sv \text{ (nmol/}\mu\text{l; or } \mu\text{mol/ml; or mM)}$$

Where:

Ay is the amount of Isocitrate (nmol) in your sample from the Standard Curve.

Sv is the sample volume (µl) added to the sample well.

Isocitrate molecular weight: 192.12 g/mol



Isocitrate Standard Curve generated using this kit protocol.

**RELATED PRODUCTS:**

- |                                     |                                      |
|-------------------------------------|--------------------------------------|
| Apoptosis Detection Kits & Reagents | Cell Proliferation & Senescence Kits |
| Glucose and Sucrose Assay Kit       | Cholesterol, LDL/HDL Assay Kits      |
| Glutathione Assay Kit               | Ethanol and Uric Acid Assay Kit      |
| NAD/NADH and NADP/NADPH Assay Kit   | Lactate Assay Kits                   |
| TAC Total Antioxidant Capacity      | Mono or Polysaccharide Assay Kits    |
| Malic Acid Assay Kit                | Glycogen/Starch Assay Kit            |

**FOR RESEARCH USE ONLY! Not to be used on humans.**

## GENERAL TROUBLESHOOTING GUIDE:

Problems	Cause	Solution
Assay not working	<ul style="list-style-type: none"> <li>• Use of ice-cold assay buffer</li> <li>• Omission of a step in the protocol</li> <li>• Plate read at incorrect wavelength</li> <li>• Use of a different 96-well plate</li> </ul>	<ul style="list-style-type: none"> <li>• Assay buffer must be at room temperature</li> <li>• Refer and follow the data sheet precisely</li> <li>• Check the wavelength in the data sheet and the filter settings of the instrument</li> <li>• Fluorescence: Black plates (clear bottoms) ; Luminescence: White plates ; Colorimeters: Clear plates</li> </ul>
Samples with erratic readings	<ul style="list-style-type: none"> <li>• Use of an incompatible sample type</li> <li>• Samples prepared in a different buffer</li> <li>• Samples were not deproteinized (if indicated in datasheet)</li> <li>• Cell/ tissue samples were not completely homogenized</li> <li>• Samples used after multiple free-thaw cycles</li> <li>• Presence of interfering substance in the sample</li> <li>• Use of old or inappropriately stored samples</li> </ul>	<ul style="list-style-type: none"> <li>• Refer data sheet for details about incompatible samples</li> <li>• Use the assay buffer provided in the kit or refer data sheet for instructions</li> <li>• Use the 10 kDa spin cut-off filter or PCA precipitation as indicated</li> <li>• Use Dounce homogenizer (increase the number of strokes); observe for lysis under microscope</li> <li>• Aliquot and freeze samples if needed to use multiple times</li> <li>• Troubleshoot if needed, deproteinize samples</li> <li>• Use fresh samples or store at correct temperatures till use</li> </ul>
Lower/ Higher readings in Samples and Standards	<ul style="list-style-type: none"> <li>• Improperly thawed components</li> <li>• Use of expired kit or improperly stored reagents</li> <li>• Allowing the reagents to sit for extended times on ice</li> <li>• Incorrect incubation times or temperatures</li> <li>• Incorrect volumes used</li> </ul>	<ul style="list-style-type: none"> <li>• Thaw all components completely and mix gently before use</li> <li>• Always check the expiry date and store the components appropriately</li> <li>• Always thaw and prepare fresh reaction mix before use</li> <li>• Refer datasheet &amp; verify correct incubation times and temperatures</li> <li>• Use calibrated pipettes and aliquot correctly</li> </ul>
Readings do not follow a linear pattern for Standard curve	<ul style="list-style-type: none"> <li>• Use of partially thawed components</li> <li>• Pipetting errors in the standard</li> <li>• Pipetting errors in the reaction mix</li> <li>• Air bubbles formed in well</li> <li>• Standard stock is at an incorrect concentration</li> <li>• Calculation errors</li> <li>• Substituting reagents from older kits/ lots</li> </ul>	<ul style="list-style-type: none"> <li>• Thaw and resuspend all components before preparing the reaction mix</li> <li>• Avoid pipetting small volumes</li> <li>• Prepare a master reaction mix whenever possible</li> <li>• Pipette gently against the wall of the tubes</li> <li>• Always refer the dilutions in the data sheet</li> <li>• Recheck calculations after referring the data sheet</li> <li>• Use fresh components from the same kit</li> </ul>
Unanticipated results	<ul style="list-style-type: none"> <li>• Measured at incorrect wavelength</li> <li>• Samples contain interfering substances</li> <li>• Use of incompatible sample type</li> <li>• Sample readings above/below the linear range</li> </ul>	<ul style="list-style-type: none"> <li>• Check the equipment and the filter setting</li> <li>• Troubleshoot if it interferes with the kit</li> <li>• Refer data sheet to check if sample is compatible with the kit or optimization is needed</li> <li>• Concentrate/ Dilute sample so as to be in the linear range</li> </ul>

**Note:** The most probable list of causes is under each problem section. Causes/ Solutions may overlap with other problems.