

Ascorbic Acid Colorimetric/Fluorometric Assay Kit

(Catalog #K661-100; 100 assays; Store at -20°C)

I. Introduction:

Ascorbic Acid (Vitamin C) plays an important role in many biological processes. It is a potent anti-oxidant, anti-inflammatory, anti-viral agent, and an immune stimulant and is present in a wide variety of foods and biological specimens. It is important to be able to monitor ascorbic acid content in these different samples. BioVision's Ascorbic Acid Assay Kit provides a rapid, simple, and sensitive means of detecting ascorbic acid in various biological samples. In this assay, our proprietary catalyst oxidizes ascorbic acid to produce a product that interacts with the ascorbic acid probe, generating color and fluorescence. Ascorbic acid can be easily determined by either colorimetric (spectrophotometry at $\lambda = 570$ nm) or fluorometric (Ex/Em = 535/587 nm) methods. The assay can detect 0.01-10 nmol of ascorbic acid per assay in various samples.

II. Kit Contents:

Components	K661-100	Cap Code	Part Number
Ascorbic Acid Assay Buffer	25 ml	WM	K661-100-1
Ascorbic Acid Probe (in DMSO)	0.2 ml	Red	K661-100-2A
Catalyst	0.5 ml	Blue	K661-100-4
Ascorbic Acid Enzyme Mix (lyophilized)	1 vial	Green	K661-100-5
Ascorbic Acid Standard (20 μ mole)	1vial	Yellow	K661-100-6

III. Storage and Handling:

Store kit at -20°C, protect from light. Warm Ascorbic Acid Assay Buffer to room temperature before use. Briefly centrifuge all small vials prior to opening.

IV. Reagent Preparation:

Ascorbic Probe: Ready to use as supplied. Warm to room temperature prior to use to completely melt frozen DMSO, then vortex to ensure uniformity. Store at -20°C, protect from light and moisture. Use within two months.

Ascorbic Acid Enzyme Mix: Dissolve in 220 μ l Ascorbic Acid Assay Buffer. Aliquot and store at -20°C. Use within two months.

Ascorbic Standard: Dissolve in 200 μ l of distilled water to generate 100 mM Ascorbic Standard stock solution. Store at -20°C. Use within two months.

Catalyst: Ready to use as supplied

V. Ascorbic Acid Assay Protocol:

1. Standard Curve Preparations:

For the colorimetric assay, dilute the standard to 1 mM by adding 10 μ l of the 100 mM Ascorbic Acid Standard to 990 μ l of distilled water, mix well. Add 0, 2, 4, 6, 8, 10 μ l into each well individually. Adjust volume to 120 μ l/well with Ascorbic Acid Assay Buffer to generate 0, 2, 4, 6, 8, 10 nmol/well of Ascorbic Acid Standard.

For the fluorometric assay, dilute the Ascorbic Acid Standard to 0.01- 0.1 mM with the Ascorbic Acid Assay Buffer (**Note:** Detection sensitivity is 10 to 100 fold higher for a fluorometric than a colorimetric assay). Follow the procedure for the colorimetric assay.

Note: Diluted ascorbic acid standard is unstable, use fresh dilution each time.

2. **Sample Preparation:** Prepare test samples to a final volume of 120 μ l/well with Ascorbic Acid Assay Buffer in a 96-well plate. We suggest testing several doses of your sample to make sure the readings are within the standard curve range.

NOTES:

1) Due to high protein content and other compounds present in serum we recommend using FRASC Ascorbic Acid Kit (K671-100) for serum samples.
2) Ascorbate is easily oxidized during sample preparation and great care must be exercised to achieve quantitative recovery.

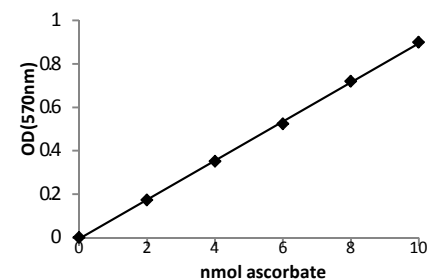
- Catalyst:** Add 100 μ l of catalyst to 900 μ l of distilled water and vortex well.
- Add 30 μ l of catalyst to each standard and sample well.
- Ascorbic Acid Reaction Mix:** Mix enough reagent for the number of samples and standards to be performed: For each well, prepare a total 50 μ l Reaction Mix containing:

46 μ l Ascorbic Acid Assay Buffer
2 μ l Ascorbic Acid Probe
2 μ l Ascorbic Acid Enzyme Mix

- Mix well. Add 50 μ l of the Reaction Mix to each well containing the Ascorbic Acid Standard and test samples. Mix well.
- Protect from light, Color is developed within 3 min and stable for an hour.
- Measure OD 570nm for colorimetric assay or Ex/Em = 535/590 nm for fluorometric assay in a micro-plate reader.
- Correct background by subtracting the value derived from the 0 ascorbic acid standard from all sample readings (Note: The background reading can be significant and must be subtracted from sample readings). Apply sample readings to the generated standard curve. Ascorbic Acid concentration can then be calculated:

$$C = As / Sv \text{ nmol}/\mu\text{l or } \mu\text{mol}/\text{ml or mM}$$

Where: **As** is ascorbic acid amount from standard curve (nmol).
Sv is the sample volume added in sample wells (μ l).
Ascorbic Acid molecular weight: 176.12.



RELATED PRODUCTS

Apoptosis Detection Kits & Reagents	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
Pyruvate and Lactate Assay Kits	cAMP/cGMP Kits

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

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