

ab65346

Ascorbic Acid Assay Kit

Instructions for Use

For the rapid, sensitive and accurate measurement of Ascorbic Acid in various samples

This product is for research use only and is not intended for diagnostic use.

Table of Contents

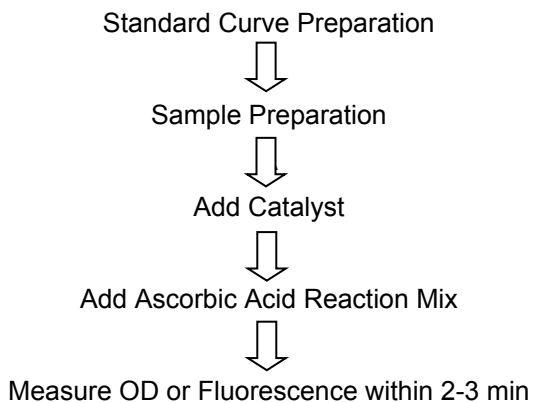
1. Overview	3
2. Protocol Summary	4
3. Components and Storage	5
4. Assay Protocol	7
5. Data Analysis	10
6. Troubleshooting	11

1. Overview

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

Abcam'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 OxiRed™ 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.

2. Protocol Summary



Components and Storage

A. Kit Components

Item	Quantity
Assay Buffer 2	25 mL
OxiRed™ Probe	0.2 mL
Catalyst Reagent	500 µL
Developer Solution V	1 vial
Ascorbic Acid Standard	1 vial

PLEASE NOTE: Assay Buffer 2 was previously labelled as Assay Buffer II and Ascorbate Acid Buffer, and Developer Solution V as Ascorbic Acid Enzyme Mix (Lyophilized), and OxiRed™ Probe as Ascorbic Acid Probe (DMSO), and Catalyst Reagent as Catalyst. The composition has not changed.

* Store kit at -20°C, protect from light. Warm Assay Buffer 2 to room temperature before use. Briefly centrifuge all small vials prior to opening.

OxiRed™ 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.

Developer Solution V: Dissolve in 220 μl Assay Buffer 2. Aliquot and store at -20°C . Use within two months.

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

Catalyst Reagent: Ready to use as supplied

B. Additional Materials Required

- Microcentrifuge
- Pipettes and pipette tips
- Fluorescent or colorimetric microplate reader
- 96 well plate
- Orbital shaker

3. Assay Protocol

1. Standard Curve Preparation:

a. For the colorimetric assay:

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

b. For the fluorometric assay:

Dilute the Ascorbic Acid Standard to 0.01- 0.1 mM with the Assay Buffer 2 (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 μl /well with Assay Buffer 2 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:

- a) Due to high protein content and other compounds present in plasma and serum we recommend using ab65656 (Ascorbic Acid Assay Kit (Biological Samples)) for plasma and serum samples.
- b) Ascorbate is easily oxidized during sample preparation and great care must be exercised to achieve quantitative recovery.

3. Add 100 μl of Catalyst Reagent to 900 μl of distilled water and vortex well. Add 30 μl of Catalyst Reagent to each standard and sample well.

4. Ascorbic Acid Reaction Mix: Mix enough reagent for the number of samples and standards to be performed: For each well, prepare a total 50 μl Reaction Mix containing:

Assay Buffer 2	46 μl
OxiRed™ Probe	2 μl
Developer Solution V	2 μl

The fluorometric assay is ~10 times more sensitive than the colorimetric assay. Use 0.4 μl of the OxiRed™ Probe per reaction to decrease the background reading/increase detection sensitivity significantly.

Mix well. Add 50 μl of the Reaction Mix to each well containing the Ascorbic Acid Standard and test samples. Mix well.

Note: Protect from light, Color is developed within 3 min and stable for an hour.

5. Measure OD_{570nm} for colorimetric assay or Ex/Em = 535/590 nm for fluorometric assay in a micro-plate reader.

Data Analysis

Correct background by subtracting the value derived from the zero ascorbic acid standard from all sample readings. 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:

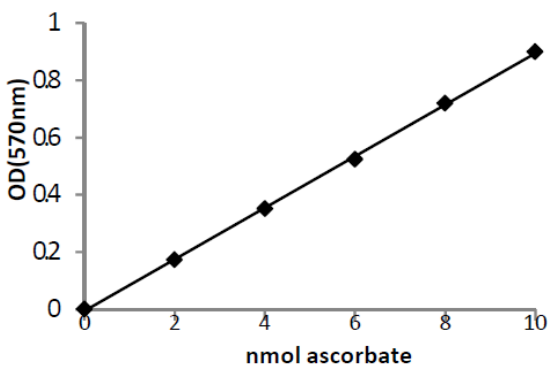
$$\text{Concentration} = \text{As} / \text{Sv} \text{ (nmol/}\mu\text{l or } \mu\text{mol/ml or mM)}$$

Where:

As is ascorbic acid amount from standard curve (nmol).

Sv is the sample volume added in sample wells (μl).

Ascorbic Acid molecular weight: 176.12.



Standard Curve performed according to assay protocol.

4. Troubleshooting

Problem	Reason	Solution
Assay not working	Assay buffer at wrong temperature	Assay buffer must not be chilled - needs to be at RT
	Protocol step missed	Re-read and follow the protocol exactly
	Plate read at incorrect wavelength	Ensure you are using appropriate reader and filter settings (refer to datasheet)
	Unsuitable microtiter plate for assay	Fluorescence: Black plates (clear bottoms); Luminescence: White plates; Colorimetry: Clear plates. If critical, datasheet will indicate whether to use flat- or U-shaped wells
Unexpected results	Measured at wrong wavelength	Use appropriate reader and filter settings described in datasheet
	Samples contain impeding substances	Troubleshoot and also consider deproteinizing samples
	Unsuitable sample type	Use recommended samples types as listed on the datasheet
	Sample readings are outside linear range	Concentrate/ dilute samples to be in linear range

Samples with inconsistent readings	Unsuitable sample type	Refer to datasheet for details about incompatible samples
	Samples prepared in the wrong buffer	Use the assay buffer provided (or refer to datasheet for instructions)
	Samples not deproteinized (if indicated on datasheet)	Use the 10kDa spin column (ab93349)
	Cell/ tissue samples not sufficiently homogenized	Increase sonication time/ number of strokes with the Dounce homogenizer
	Too many freeze-thaw cycles	Aliquot samples to reduce the number of freeze-thaw cycles
	Samples contain impeding substances	Troubleshoot and also consider deproteinizing samples
	Samples are too old or incorrectly stored	Use freshly made samples and store at recommended temperature until use
Lower/ Higher readings in samples and standards	Not fully thawed kit components	Wait for components to thaw completely and gently mix prior use
	Out-of-date kit or incorrectly stored reagents	Always check expiry date and store kit components as recommended on the datasheet
	Reagents sitting for extended periods on ice	Try to prepare a fresh reaction mix prior to each use
	Incorrect incubation time/ temperature	Refer to datasheet for recommended incubation time and/ or temperature
	Incorrect amounts used	Check pipette is calibrated correctly (always use smallest volume pipette that can pipette entire volume)

Problem	Reason	Solution
Standard curve is not linear	Not fully thawed kit components	Wait for components to thaw completely and gently mix prior use
	Pipetting errors when setting up the standard curve	Try not to pipette too small volumes
	Incorrect pipetting when preparing the reaction mix	Always prepare a master mix
	Air bubbles in wells	Air bubbles will interfere with readings; try to avoid producing air bubbles and always remove bubbles prior to reading plates
	Concentration of standard stock incorrect	Recheck datasheet for recommended concentrations of standard stocks
	Errors in standard curve calculations	Refer to datasheet and re-check the calculations
	Use of other reagents than those provided with the kit	Use fresh components from the same kit

For further technical questions please do not hesitate to contact us by email (technical@abcam.com) or phone (select “*contact us*” on www.abcam.com for the phone number for your region).

Technical Support

Copyright © 2025 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

www.abcam.com/contactus

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)