

ab155900

Xanthine/Hypoxanthine Assay Kit

Instructions for Use

For the sensitive and accurate measurement of Xanthine/Hypoxanthine in various tissues/cells and body fluids and the analysis of purine metabolism and cell signaling.

[View kit datasheet: www.abcam.com/ab155900](http://www.abcam.com/ab155900)

(use www.abcam.cn/ab155900 for China, or www.abcam.co.jp/ab155900 for Japan)

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

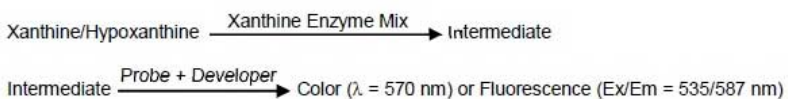
PLEASE NOTE: With the acquisition of BioVision by Abcam, we have made some changes to component names and packaging to better align with our global standards as we work towards environmental-friendly and efficient growth. You are receiving the same high-quality products as always, with no changes to specifications or protocols.

Table of Contents

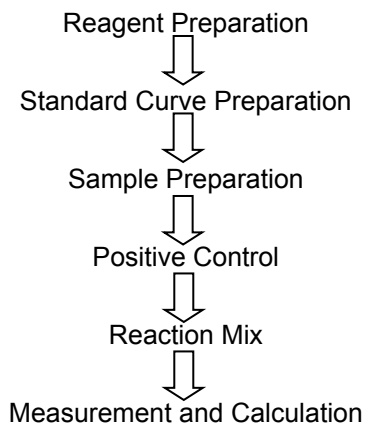
| | |
|---------------------------|----|
| 1. Overview | 3 |
| 2. Protocol Summary | 4 |
| 3. Components and Storage | 5 |
| 4. Assay Protocol | 7 |
| 5. Data Analysis | 10 |
| 6. Troubleshooting | 12 |

1. Overview

Xanthine, a catabolic product of purine metabolism, is present in body fluids, muscle tissue and certain plants. Structurally like caffeine, Xanthine has a stimulant effect and is used clinically to treat the congestive diseases such as asthma and chronic obstructive pulmonary disease. Xanthine is metabolized into uric acid and superoxide by Xanthine oxidase. Xanthine oxidase deficiency causes the rare genetic disorder-Xanthinuria, and leads to Xanthine accumulation in urine and blood, which ultimately progresses to renal failure. Recent studies show that Xanthine levels are elevated following ischemic injury, thus Xanthine can serve as a useful marker for tissue hypoxia. Early detection of Xanthine alteration in biological fluids is crucial for metabolic studies and for diagnostic and therapeutic monitoring. In Abcam's Xanthine/Hypoxanthine Assay kit, Xanthine/Hypoxanthine is specifically oxidized by the Xanthine Development Enzyme Mix VIII/Enzyme Mix to form an intermediate, which reacts with Developer Solution V/Developer & Probe to form a product that can be measured colorimetrically ($\lambda = 570 \text{ nm}$) or fluorometrically (Ex/Em = 535/587 nm). Xanthine/Hypoxanthine Assay kit is rapid, simple and sensitive. This high-throughput suitable assay kit can detect Xanthine levels as low as $0.4 \mu\text{M}$ in various biological samples.



2. Protocol Summary



3. Components and Storage

A. Kit Components

| Item | Quantity |
|--|----------|
| Assay Buffer II/Assay Buffer | 25 mL |
| OxiRed Probe/OxiRed™ Probe (in DMSO) | 0.2 mL |
| Development Enzyme Mix VIII/Enzyme Mix (Lyophilized) | 1 vial |
| Developer Solution V/Developer (Lyophilized) | 1 vial |
| Xanthine Standard/Xanthine Standard (Lyophilized) | 1 vial |

* Store the kit at -20°C and protect from light. Please read the entire protocol before performing the assay. Avoid repeated freeze/thaw cycles.

Warm all Buffers to room temperature before use. Briefly centrifuge all small vials prior to opening.

B. Additional Materials Required

- 96-well clear plate with flat bottoms (colorimetric)
- 96-well white plate with flat bottoms (fluorometric)
- Multi-well spectrophotometer (ELISA reader)

4. Assay Protocol

A. Reagent Preparation

1. Development Enzyme Mix VIII/Enzyme Mix:

Reconstitute with 220 μ l Assay Buffer II/Assay Buffer. Pipette up and down to dissolve completely. Aliquot and store at -20°C . Avoid repeated freeze/thaw cycles. Use within two months. Keep on ice while in use.

2. Developer Solution V/Developer:

Reconstitute with 220 μ l Assay Buffer II/Assay Buffer. Pipette up and down to dissolve completely. Aliquot and store at -20°C . Use within two months. Keep on ice while in use.

3. Xanthine Standard:

Reconstitute with 500 μ l dH_2O to generate 2.0 mM (2.0 nmol/ μ l) Xanthine Standard solution. Store at -20°C . Use within two months. Keep on ice while in use.

B. Xanthine Assay Protocol

1. Xanthine Standard Curve:

For colorimetric assay, add 0, 2, 4, 6, 8 & 10 μ l of 2 mM Xanthine Standard into series of wells in 96 well plate to generate 0, 4, 8, 12, 16 & 20 nmol/well Xanthine Standard.

Adjust volume to 50 μ l per well with Assay Buffer II/Xanthine Assay Buffer.

For fluorometric assay, dilute Xanthine Standard to 0.02 mM (20 pmol/ μ l) by adding 10 μ l of 2 mM Xanthine Standard to 990 μ l dH₂O & mix. Add 0, 2, 4, 6, 8 & 10 μ l of 0.02 mM Xanthine Standard into series of wells in 96 well plate to generate 0, 40, 80, 120, 160 & 200 pmol/well Xanthine Standard. Adjust volume to 50 μ l per well with Assay Buffer II/Xanthine Assay Buffer.

2. Sample Preparation:

Liquid samples can be measured directly. Rapidly homogenize tissue (10 mg) or cells (1×10^6) with 100 μ l ice cold Assay Buffer II/Assay Buffer for 10 minutes on ice. Centrifuge at 12000 rpm for 5 min. Collect the supernatant. Add 1-50 μ l sample per well, adjust final volume to 50 μ l with Assay Buffer II/Assay Buffer.

Notes:

- a) *Some enzymes in samples may interfere with the assay. Enzymes can be removed by 10 K quick spin columns or by perchloric acid/KOH treatment.*
- b) *For unknown samples, we suggest testing several doses to ensure the readings are within the Standard Curve range.*
- c) *For samples having high background, prepare parallel sample well(s) as background control.*

3. Reaction Mix:

Mix enough reagents for the number of assays to be performed.

For each well, prepare 50 μ l Reaction Mix containing:

| | Reaction Mix | Background Control Mix |
|--------------------|---------------------|-------------------------------|
| Assay Buffer | 44 μ l | 46 μ l |
| II/Assay Buffer | | |
| Development | 2 μ l | --- μ l |
| Enzyme Mix | | |
| VIII/Enzyme Mix | | |
| Developer Solution | 2 μ l | 2 μ l |
| V/Developer | | |
| Probe | 2 μ l | 2 μ l |

Add 50 μ l of the Reaction Mix to each well containing the Standard and test. Mix well.

Note: For samples having high background, add 50 μ l of the Background Control Mix to sample background control well(s). Mix well

4. Measurement:

Incubate for 30 min at room temperature, protected from light. Measure fluorescence at Ex/Em = 535/587 nm or color at $\lambda = 570$ nm.

5. Data Analysis

Calculation: Subtract 0 Standard reading from all readings. Plot the Xanthine Standard Curve. For samples having high background, correct sample background by subtracting the value derived from the background control from sample readings. Apply the corrected sample reading to the Xanthine Standard Curve to get B pmol or nmol of Xanthine/Hypoxanthine in the sample(s).

$$\text{Xanthine/Hypoxanthine Concentration in sample} = \frac{\mathbf{B}}{\mathbf{V}} \times \text{Dilution Factor} = \text{nmol/ml/ or pmol/ml} = \mu\text{M or nM}$$

Where:

B is the amount of Xanthine/Hypoxanthine in the sample (pmol or nmol).

V is the sample volume added into the reaction well (ml).

Xanthine molecular weight: 152.11 g/mol. Hypoxanthine molecular weight: 136.11g/mol

Xanthine/Hypoxanthine in samples can also be expressed in nmol/mg of sample or other desired method.

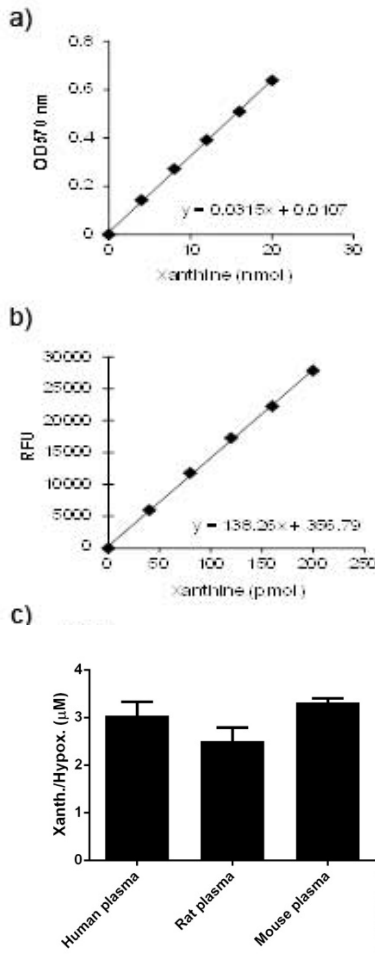


Figure 1: Xanthine Standard Curve (a) & (b). Xanthine & Hypoxanthine measured in biological fluids by fluorometric method showing concentrations (micromolar) (c). Assays were performed following Kit protocol.

6. 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 |

| Problem | Reason | Solution |
|---|---|---|
| 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 © 2023 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)