

# Xanthine Oxidase Activity Assay Kit ab102522

[19 References](#) [2 Images](#)

### Overview

<b>Product name</b>	Xanthine Oxidase Activity Assay Kit
<b>Detection method</b>	Colorimetric/Fluorometric
<b>Sample type</b>	Milk, Urine, Serum, Plasma, Cell culture extracts, Other biological fluids, Tissue Extracts
<b>Assay type</b>	Enzyme activity
<b>Sensitivity</b>	> 10 mU/ml
<b>Range</b>	10 mU/ml - 1000 mU/ml
<b>Assay time</b>	0h 40m
<b>Product overview</b>	<p>Xanthine Oxidase Activity Assay Kit ab102522 is an easy and sensitive assay to determine xanthine oxidase activity in variety of samples.</p> <p>In the xanthine oxidase assay protocol, xanthine oxidase oxidizes xanthine to hydrogen peroxide (<math>H_2O_2</math>) which reacts stoichiometrically with a probe to generate color (at OD = 570 nm) and fluorescence (at Ex/Em = 535/587 nm). Since the color or fluorescence intensity is proportional to XO content, the XO activity can be accurately measured.</p> <p>The kit detects 1-100 mU xanthine oxidase in 100 <math>\mu</math>l reaction volume.</p> <p>Xanthine oxidase assay protocol summary:</p> <ul style="list-style-type: none"> <li>- add samples and standards to wells</li> <li>- add reaction mix</li> <li>- analyze with microplate reader immediately and after 10-20 min</li> </ul>
<b>Notes</b>	<p>This product is manufactured by BioVision, an Abcam company and was previously called K710 Xanthine Oxidase Activity Colorimetric/Fluorometric Assay Kit. K710-100 is the same size as the 100 test size of ab102522.</p> <p>Xanthine oxidase (XO, EC 1.17.3.2 ) is present in appreciable amounts in the liver and jejunum in healthy individuals. However, in various liver disorders, XO is released into circulation. Therefore, determination of serum XO level serves as a sensitive indicator of acute liver damage such as jaundice.</p>
<b>Platform</b>	Microplate reader

### Properties

**Storage instructions**

Store at -20°C. Please refer to protocols.

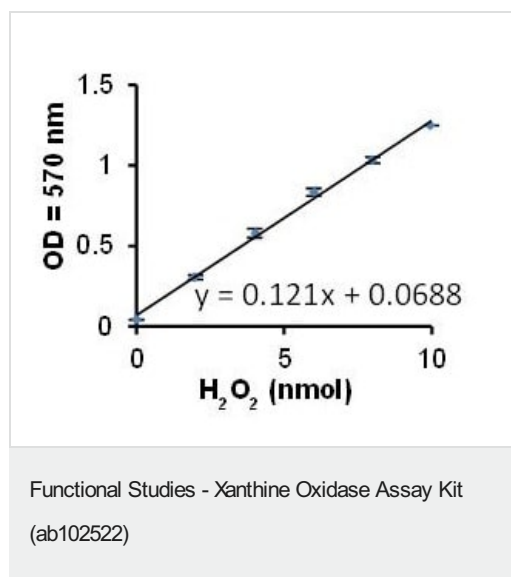
Components	100 tests
Assay Buffer VII	1 x 25ml
Developer Solution V	1 vial
H <sub>2</sub> O <sub>2</sub> Standard	1 x 0.1ml
OxiRed Probe	1 x 200µl
Xanthine Enzyme Mix	1 vial
XO Substrate Mix	1 vial

**Relevance**

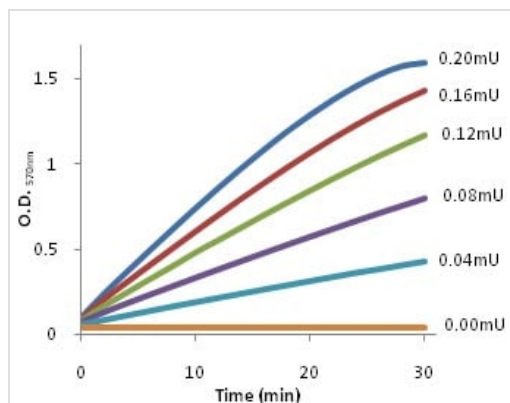
Xanthine oxidase (XO, EC 1.17.3.2 ) is present in appreciable amounts in the liver and jejunum in healthy individuals. However, in various liver disorders, XO is released into circulation. Therefore, determination of serum XO level serves as a sensitive indicator of acute liver damage such as jaundice.

**Cellular localization**

Cytoplasm. Peroxisome. Secreted.

**Images**

Representative Standard Curve using ab102522.



Xanthine Oxidase Positive Control.

Functional Studies - Xanthine Oxidase Assay Kit  
(ab102522)

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

### Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

### Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors