

Peroxidase Activity Assay Kit ab155895

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

Overview

Product name	Peroxidase Activity Assay Kit
Detection method	Colorimetric/Fluorometric
Sample type	Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids
Assay type	Enzyme activity
Species reactivity	Reacts with: Mammals, Other species
Product overview	<p>Abcam's Peroxidase Assay Kit (ab155895) provides a convenient colorimetric and fluorometric means to measure the peroxidase activity in biological samples. In the presence of Peroxidase, the OxiRed Probe reacts with H₂O₂ in a 1:1 stoichiometry to produce the redfluorescent oxidation product, resorufin. The resorufin is quantified by colorimetric ($\lambda_{\text{max}} = 570\text{nm}$) or fluorometric methods ($\text{Ex/Em} = 535/587\text{ nm}$). The assay is simple, direct, highly sensitive and high throughput-ready. The detection limit is 0.1 mU per well via colorimetric or 0.01 mU per well via fluorometric method, based on our unit definition.</p>
Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K772 Peroxidase Activity Colorimetric/Fluorometric Assay Kit. K772-100 is the same size as the 100 test size of ab155895.</p> <p>Peroxidases (EC number 1.11.1.x) are a large family of enzymes that typically catalyze a reaction of the form: ROOR' + electron donor (2 e-) + 2H⁺ → ROH + R'OH. For many of these enzymes the optimal substrate is hydrogen peroxide, but others are more active with organic hydroperoxides such as lipid peroxides. Peroxidases can contain a heme cofactor in their active sites, or alternately redox-active cysteine or selenocysteine residues.</p>
Platform	Microplate reader

Properties

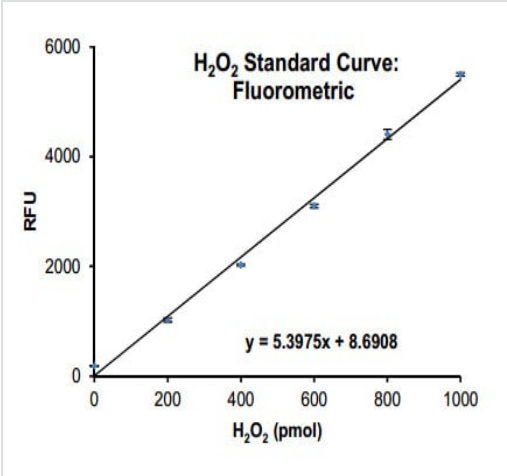
Storage instructions Store at -20°C. Please refer to protocols.

Components	100 tests
Assay Buffer XVIII	1 x 25ml
Developer Solution V	1 vial

Components	100 tests
H2O2 Standard	1 x 0.1ml
OxiRed Probe	1 x 0.2ml

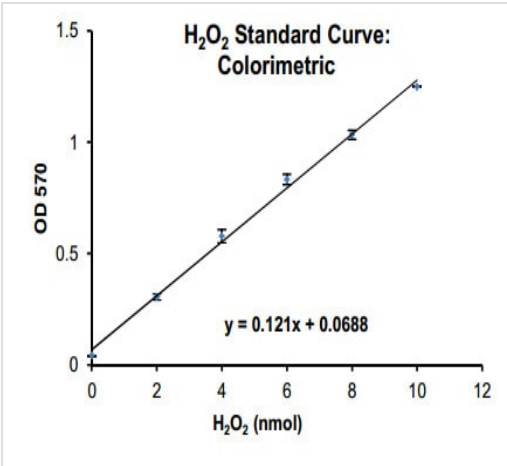
Relevance	<p>Horseradish Peroxidase (HRP) is an enzyme commonly used as an indicator for chemical reactions which produce peroxide. The enzyme is routinely conjugated to antibodies for use in enzyme-based immunoassay systems. HRP functions in the removal of H2O2 (hydrogen peroxide), oxidation of toxic reductants, biosynthesis and degradation of lignin, suberization, auxin catabolism, response to environmental stresses such as wounding, pathogen attack and oxidative stress. These functions might be dependent on each isozyme/isoform in each plant tissue.</p>
Cellular localization	<p>Secreted Probable. Vacuole Probable. Note: Carboxy-terminal extension appears to target the protein to vacuoles.</p>

Images



An example of a Fluorimetric Standard Curve

H2O2 standard curve - fluorometric



An example of a Colorimetric Standard Curve

H2O2 Standard Curve - Colorimetric

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