

Enolase Assay Kit ab241024

[1 References](#) [3 Images](#)

Overview

Product name	Enolase Assay Kit
Detection method	Colorimetric/Fluorometric
Sample type	Cell culture supernatant, Tissue, Adherent cells, Suspension cells
Product overview	In the Enolase Assay Kit (ab241024) enolase catalyzes the conversion of 2-phosphoglycerate to phosphoenolpyruvate, which is subsequently used to generate an intermediate product. The intermediate product stoichiometrically reacts with the probe to generate color (OD 570 nm) or fluorescence (Ex/Em = 535/587 nm).

This simple & sensitive assay Kit can detect enolase activity less than 0.04 mU in a variety of samples.

Notes	This product is manufactured by BioVision, an Abcam company and was previously called K691 Enolase Activity Colorimetric/Fluorometric Assay Kit. K691-100 is the same size as the 100 test size of ab241024.
--------------	--

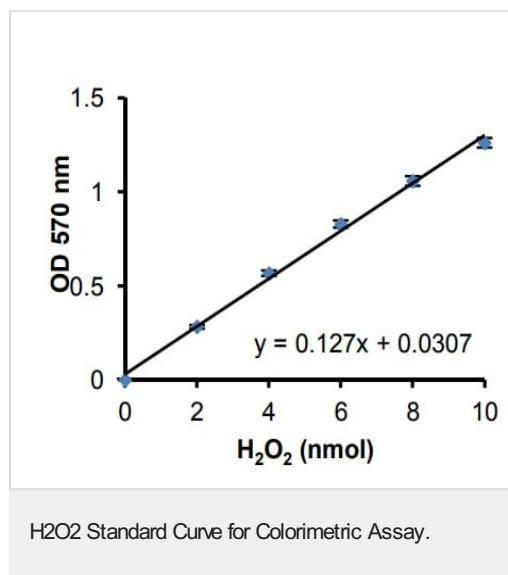
Properties

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

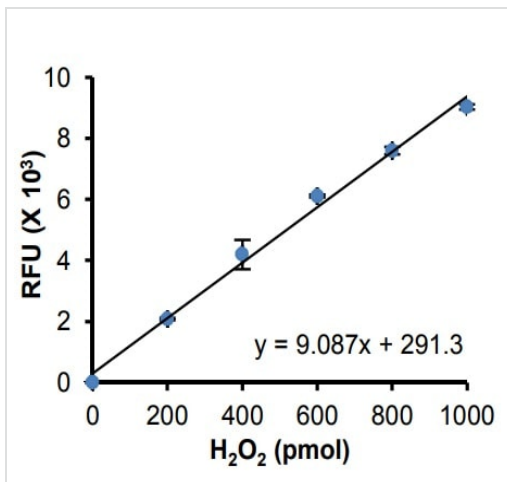
Components	100 tests
2-Phosphoglycerate	1 vial
Assay Buffer IV	1 x 25ml
Developer V	1 vial
Enolase Converter Mix	1 vial
Enolase Positive Control	1 vial
H2O2 Standard	1 x 0.1ml
OxiRed Probe	1 x 0.2ml

Function	Multifunctional enzyme that, as well as its role in glycolysis, plays a part in various processes such as growth control, hypoxia tolerance and allergic responses. May also function in the intravascular and pericellular fibrinolytic system due to its ability to serve as a receptor and activator of plasminogen on the cell surface of several cell-types such as leukocytes and neurons. Stimulates immunoglobulin production. MBP1 binds to the myc promoter and acts as a transcriptional repressor. May be a tumor suppressor.
Tissue specificity	The alpha/alpha homodimer is expressed in embryo and in most adult tissues. The alpha/beta heterodimer and the beta/beta homodimer are found in striated muscle, and the alpha/gamma heterodimer and the gamma/gamma homodimer in neurons.
Pathway	Carbohydrate degradation; glycolysis; pyruvate from D-glyceraldehyde 3-phosphate: step 4/5.
Sequence similarities	Belongs to the enolase family.
Developmental stage	During ontogenesis, there is a transition from the alpha/alpha homodimer to the alpha/beta heterodimer in striated muscle cells, and to the alpha/gamma heterodimer in nerve cells.
Post-translational modifications	ISGylated.
Cellular localization	Nucleus and Cytoplasm. Cell membrane. Cytoplasm > myofibril > sarcomere > M line. Can translocate to the plasma membrane in either the homodimeric (alpha/alpha) or heterodimeric (alpha/gamma) form. ENO1 is localized to the M line.

Images

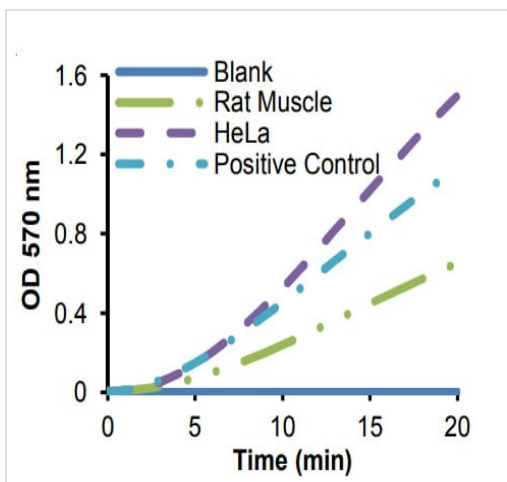


H₂O₂ Standard Curve for Colorimetric Assay.



H₂O₂ Standard Curve for Fluorometric Assay.

H₂O₂ Standard Curve for Fluorometric Assay.



Enolase Activity in rat muscle lysate (1 µg), HeLa lysate (5 µg) and positive control. Assays were performed following the kit protocol.

Enolase Activity.

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