

Alpha-Glucosidase Activity Assay Kit (Colorimetric) ab174093

2 References [画像数 4](#)

製品の概要

| | |
|---------|--|
| 製品名 | Alpha-Glucosidase Activity Assay Kit (Colorimetric) |
| 検出方法 | Colorimetric |
| サンプルの種類 | Saliva, Serum, Cell culture extracts, Tissue, Adherent cells, Suspension cells |
| アッセイタイプ | Enzyme activity (quantitative) |
| 検出範囲 | 0.1 mU/well - 10 mU/well |
| 種交差性 | 交差種: Mammals, Other species |
| 製品の概要 | In Abcam's Alpha-Glucosidase Activity Assay Kit (Colorimetric) (ab174093), α -Glucosidase hydrolyzes the Substrate Mix to release the p-nitrophenol that can be measured colorimetrically (OD = 410 nm). This is an easy, quick and high-throughput capable kit that can measure 0.1-10 mU of α -glucosidase activity in a variety of samples. |

Visit our [FAQs page](#) for tips and troubleshooting.

| | |
|------|--|
| 特記事項 | <p>This product is manufactured by BioVision, an Abcam company and was previously called K690 α-Glucosidase Activity Colorimetric Assay Kit. K690-100 is the same size as the 100 test size of ab174093.</p> <p>α-Glucosidase breaks down α-1,4 linked polysaccharides to glucose, which can be utilized as a source of energy. In the biotechnology industry, α-glucosidase is used to produce glucose from intermediate breakdown products of starch hydrolysis generated by enzymes such as amylase.</p> <p>Pompe disease, one of the 12 known glycogen storage diseases, is an autosomal recessive metabolic disorder attributed to α-glucosidase deficiency. In this disease, glycogen accumulates in the lysosomes, resulting in progressive muscle weakness, heart failure and other neurological symptoms.</p> |
|------|--|

| | |
|------------|-------------------|
| 試験プラットフォーム | Microplate reader |
|------------|-------------------|

製品の特性

| | |
|------|--|
| 保存方法 | Store at -20°C. Please refer to protocols. |
|------|--|

| 内容 | 100 tests |
|--------------------------------|-----------|
| α-Glucosidase Assay Buffer | 1 x 25ml |
| α-Glucosidase Positive Control | 1 vial |
| α-Glucosidase Substrate Mix | 1 x 0.3ml |
| p-Nitrophenol Standard | 1 x 100μl |

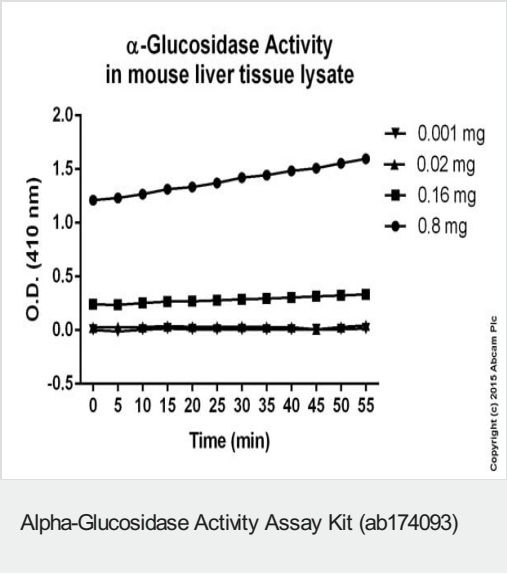
関連性

a-Glucosidase breaks down a-1,4 linked polysaccharides to glucose, which can be utilized as a source of energy. In the biotechnology industry, a-glucosidase is used to produce glucose from intermediate breakdown products of starch hydrolysis generated by enzymes such as amylase. Pompe disease, one of the 12 known glycogen storage diseases, is an autosomal recessive metabolic disorder attributed to a- glucosidase deficiency. In this disease, glycogen accumulates in the lysosomes, resulting in progressive muscle weakness, heart failure and other neurological symptoms.

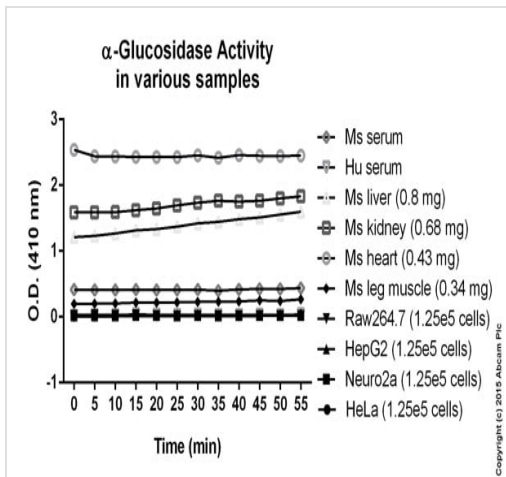
細胞内局在

Lysosome. Lysosome membrane.

画像

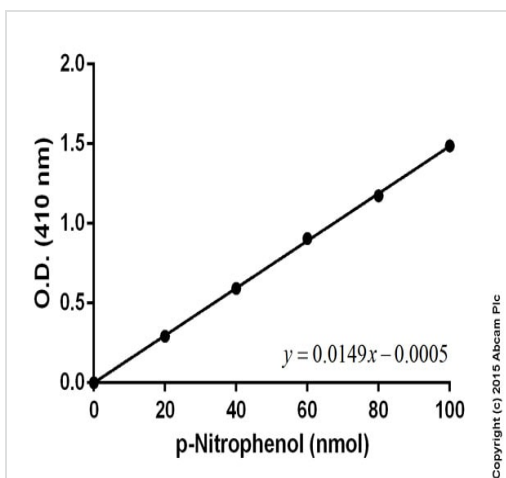


Time course of Alpha-Glucosidase Activity in mouse liver tissue lysate



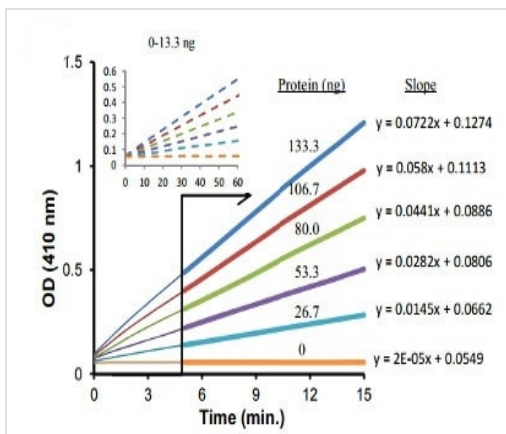
Alpha-Glucosidase Activity Assay Kit (ab174093)

Time course of Alpha-Glucosidase Activity in various samples



Alpha-Glucosidase Activity Assay Kit (ab174093)

Standard curve: mean of duplicates (+/- SD) with background reads subtracted



α-Glucosidase kinetic assay

Kinetic profile of various amounts (0, 2, 4, 6, 8 & 10 mU) of α-glucosidase run at 25°C under this protocol. Inset: Results for 0-0.2-0.4-0.6-0.8-1.0 mU of α-glucosidase. Data points after 5 minutes were used to determine slope. This is example data only.

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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