

# Aldolase Activity Assay Kit (Colorimetric) ab196994

**13 References**   **3 图像**

### 概述

|       |  |
|-------|--|
| 产品名称  | Aldolase Activity Assay试剂盒(Colorimetric)   |
| 检测方法  | Colorimetric   |
| 样品类型  | Serum, Plasma, Tissue, Adherent cells, Suspension cells  |
| 检测类型  | Enzyme activity  |
| 灵敏度   | < 100 µU   |
| 种属反应性 | 与反应: Mammals, Other species  |
| 产品概述  | Aldolase Activity Assay (Colorimetric) (ab196994), aldolase converts fructose-1,6-bisphosphate to glyceraldehyde-3-phosphate and dihydroxyacetone, and through a series of reactions, reduces a nearly colorless probe to a colored product with absorbance at 450 nm. |

This assay kit is simple, sensitive and high-throughput adaptable and it can be used to measure aldolase activity in various cells and tissues as well as to analyze glycolysis and glucogenesis pathways.

**说明** This product is manufactured by BioVision, an Abcam company and was previously called K665 Aldolase Activity Colorimetric Assay Kit. K665-100 is the same size as the 100 test size of ab196994.

Aldolase (Fructose-Bisphosphate Aldolase: EC 4.1.2.13) is an important enzyme for both glycolysis and gluconeogenesis. It catalyzes the reversible reaction of fructose-1,6-bisphosphate to glyceraldehyde-3-phosphate & dihydroxyacetone.

There are 2 classes of Aldolase - class I: found in animal and plant tissues and class II: found in prokaryotes and lower eukaryotes. Class I Aldolase has 3 isozymes- Type A: found in muscle and red blood cells, Type B: found in liver and kidney and Type C: found in brain. Aldolase A deficiency leads to myopathy & hemolytic anemia. Muscle disease and liver injury can also cause increased serum aldolase.

Accurate detection of aldolase activity is valuable for diagnostic and mechanistic studies.

**平台** Microplate reader

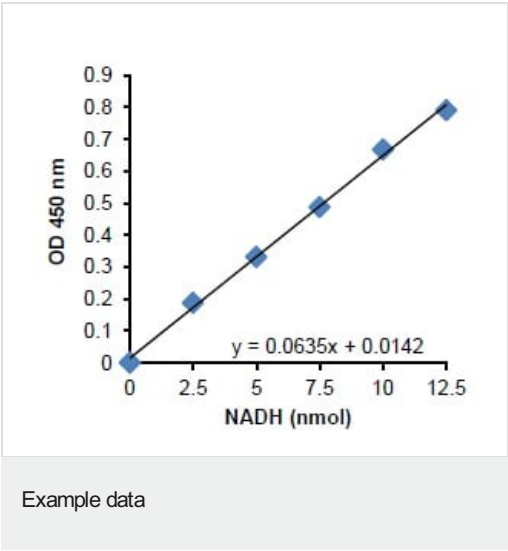
### 性能

**存放说明** Store at -20°C. Please refer to protocols.

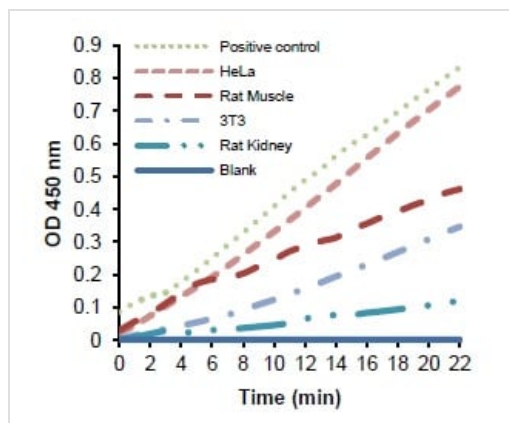
| 组件                        | 100 tests |
|---------------------------|-----------|
| PGK Developer             | 1 vial    |
| Aldolase Substrate        | 1 vial    |
| Assay Buffer IV           | 1 x 25ml  |
| Developer Solution III    | 1 vial    |
| NADH Standard I           | 1 vial    |
| Aldolase Positive Control | 1 vial    |

|       |  |
|-------|--|
| 功能    | Plays a key role in glycolysis and gluconeogenesis. In addition, may also function as scaffolding protein.   |
| 通路    | Carbohydrate degradation; glycolysis; D-glyceraldehyde 3-phosphate and glycerone phosphate from D-glucose: step 4/4.   |
| 疾病相关  | Defects in ALDOA are the cause of glycogen storage disease type 12 (GSD12) [MIM:611881]; also known as red cell aldolase deficiency. A metabolic disorder associated with increased hepatic glycogen and hemolytic anemia. It may lead to myopathy with exercise intolerance and rhabdomyolysis. |
| 序列相似性 | Belongs to the class I fructose-bisphosphate aldolase family.  |

图片

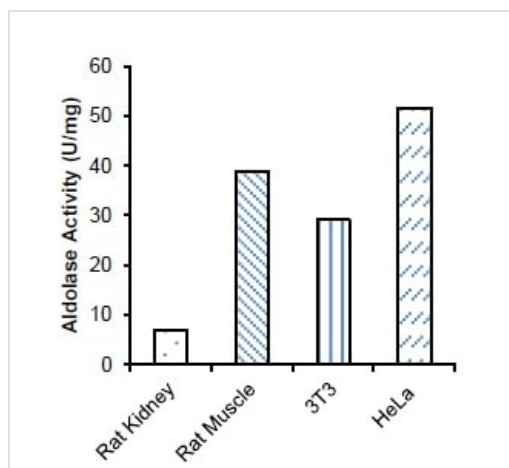


Typical NADH Standard Curve.



Example data

Aldolase activity in rat kidney and muscle, HeLa and 3T3 cells. Data obtained performing assay as described in assay protocol.



Example data

Relative aldolase activity in lysates prepared from rat kidney (14.2 µg), rat muscle (10.64 µg), 3T3 cells (9.97 µg) and HeLa cells (11.35 µg).

**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.cn/abpromise> or contact our technical team.

## Terms and conditions

---

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