

Product datasheet

Citrate Synthase Assay Kit ab239712

1 References 2 Images

Overview

<b>Product name</b>	Citrate Synthase Assay Kit
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Cell culture supernatant, Tissue, Cell culture media
<b>Product overview</b>	Citrate Synthase Assay Kit (ab239712) provides reagents for the measurement of citrate synthase activity in various tissues/cells. Citrate Synthase reacts with substrate mix to form an intermediate, which subsequently reacts with developer to generate the colored product. The rate of color development is proportional to the enzyme activity. The assay is simple, rapid and can detect Citrate Synthase activity less than 1 mU in a variety of samples.
<b>Platform</b>	Microplate reader

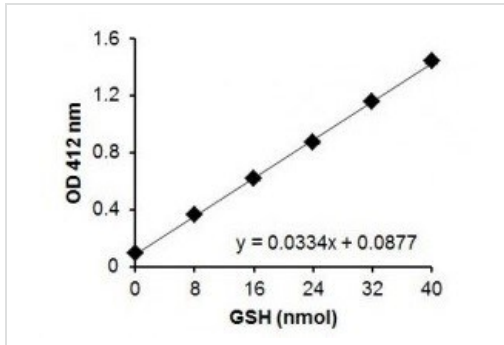
Properties

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

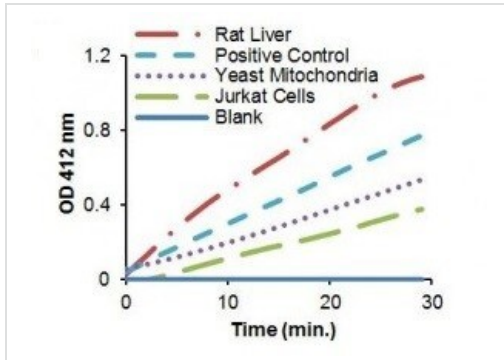
Components	100 tests
CS Assay Buffer	1 x 25ml
CS Developer (Lyophilized)	1 vial
CS Positive Control (Lyophilized)	1 vial
CS Substrate Mix (Lyophilized)	1 vial
GSH Standard (reduced) (Lyophilized)	1 vial

<b>Pathway</b>	Carbohydrate metabolism; tricarboxylic acid cycle; isocitrate from oxaloacetate: step 1/2.
<b>Sequence similarities</b>	Belongs to the citrate synthase family.
<b>Cellular localization</b>	Mitochondrion matrix.

Images



GSH Standard Curve



Citrate Synthase Activity in Jurkat cell lysate (10 µg), rat liver lysate (20 µg), purified yeast mitochondria (4 µg) and CS Positive Control (2 µL).

**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