

# Sucrose Assay Kit ab83387

[1 References](#) [1 图像](#)

### 概述

产品名称	Sucrose Assay试剂盒
检测方法	Colorimetric/Fluorometric
样品类型	Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids, Tissue Extracts
检测类型	Quantitative
灵敏度	> 0.2 $\mu$ M
范围	0 $\mu$ M - 10 $\mu$ M
检测时间	1h 00m
产品概述	Abcam's Sucrose Assay Kit provides a convenient means for specifically measuring sucrose levels from various biological samples (e.g. serum, plasma, body fluids, food, growth medium, etc.). Sucrose is converted to glucose and fructose by invertase. The generated glucose is then specifically oxidized to generate a product which reacts with a probe to generate color ( $\lambda$ = 570 nm) and fluorescence (Ex/Em = 535/587 nm). Use of proper controls allows for correcting free glucose background. Other disaccharides such as lactose and maltose do not interfere with the assay. The method can detect 0.0002-10 mM sucrose concentrations.

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

**说明** This product is manufactured by BioVision, an Abcam company and was previously called K626 Sucrose Colorimetric/Fluorometric Assay Kit. K626-100 is the same size as the 100 test size of ab83387.

Sucrose ( $C_{12}H_{22}O_{11}$ ; FW:342.3), also called table sugar, is one of the most important fuel sources to generate the universal energy molecule ATP. Sucrose is a disaccharide which can be converted into one glucose and one fructose.

**平台** Microplate reader

### 性能

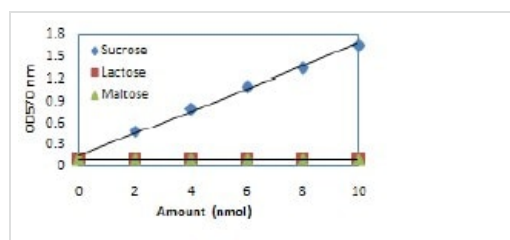
**存放说明** Store at -20°C. Please refer to protocols.

组件	100 tests
Assay Buffer II	1 x 25ml

组件	100 tests
Development Enzyme Mix II	1 vial
Invertase	1 vial
OxiRed Probe	1 x 200µl
Sucrose Standard	1 x 100µl

**相关性** Sucrose ( $C_{12}H_{22}O_{11}$ ; FW:342.3), also called table sugar, is one of the most important fuel sources to generate the universal energy molecule ATP. Sucrose is a disaccharide which can be converted into one glucose and one fructose.

## 图片



Standard Sucrose Assay. Different amounts of sucrose, lactose and maltose are assayed following the kit protocol. Data show that the Sucrose assay kit specifically detects sucrose, but not lactose or maltose.

Functional Studies - Sucrose Assay Kit (ab83387)

**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