

D-Sorbitol Assay Kit (Colorimetric) ab118968

[7 References](#) [3 图像](#)

概述

产品名称	D-Sorbitol Assay试剂盒(Colorimetric)
检测方法	Colorimetric
样品类型	Food samples
检测类型	Quantitative
灵敏度	> 0.1 nmol/well
范围	0.1 nmol/well - 10 nmol/well
检测时间	0h 40m
产品概述	Abcam's D-Sorbitol Assay Kit (Colorimetric) is designed to measure sorbitol in a variety of samples such as foods, fruits, fruit juices, pharmaceuticals, cosmetics, paper. This assay is not recommended for plasma, serum or urine samples. In the assay, sorbitol is oxidized to fructose with the proportional development of intense color with an absorbance maximum at 560 nm. The assay is useful over the range of 0.1-10 nmol of Sorbitol per sample.

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

说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K631 D-Sorbitol Colorimetric Assay Kit. K631-100 is the same size as the 100 test size of ab118968.</p> <p>Sorbitol is one of the 6 carbon sugar alcohols. It is commonly used as an artificial sweetener, as a laxative and in cosmetics as a humectant and thickening agent. Sorbitol is produced naturally in a variety of fruits. It can be produced in humans in small amounts by the reduction of glucose by aldose reductase. Due to its poor ability to diffuse across the cell membrane, sorbitol can be trapped in cells and is believed to be one of the causes of damage (due to osmotic effects) in diabetes. Interestingly, sorbitol can be used as a screen for the O154:H7 strain of E. coli, since this strain is one of the few strains which cannot metabolize sorbitol.</p>
----	--

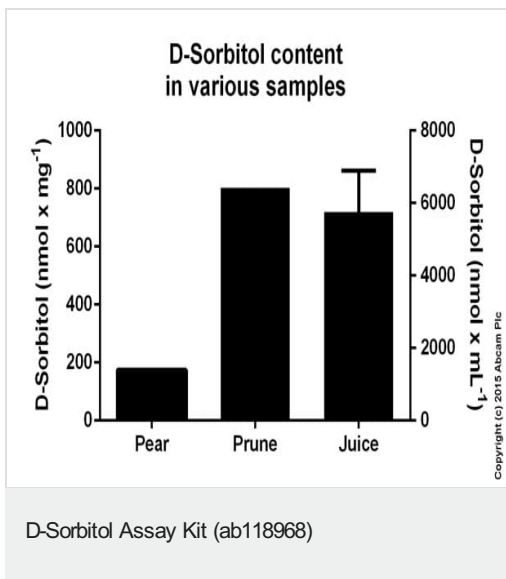
平台	Microplate reader
----	-------------------

性能

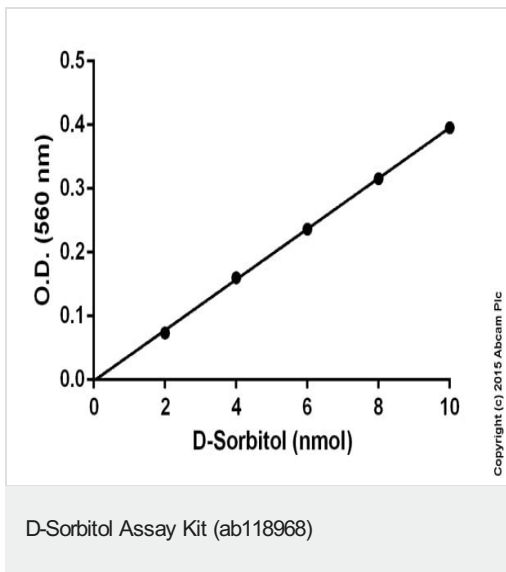
存放说明	Store at -20°C. Please refer to protocols.
------	--

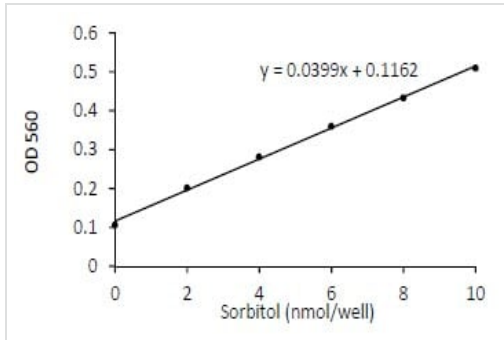
组件	100 tests
Sorbitol Assay Buffer	1 x 25ml
Sorbitol Probe	1 x 200µl
Sorbitol Developer	1 vial
Sorbitol Enzyme Mix	1 vial
Sorbitol Standard	1 x 100µl

图片



D-Sorbitol measured in various samples showing nmol/mg (fruits) and nmol/mL (juice). 300 mg (wet weight) of prune and pear was homogenised in 1 mL of water. Juice was diluted in water. Samples were diluted 2-1000 fold.





Example of standard curve obtained using ab118968.

Functional Studies - D-Sorbitol Colorimetric Assay
Kit (ab118968)

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