

## Product datasheet

# High Sensitivity Carbohydrate Assay Kit (Colorimetric) ab284543

[3 Images](#)

### Overview

---

<b>Product name</b>	High Sensitivity Carbohydrate Assay Kit (Colorimetric)
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Other biological fluids, Tissue, Adherent cells, Suspension cells, Food samples
<b>Assay type</b>	Quantitative
<b>Range</b>	0 µg - 20 µg
<b>Assay duration</b>	Multiple steps standard assay
<b>Product overview</b>	<p>High Sensitivity Carbohydrate Assay Kit (Colorimetric) (ab284543) (K2049) is a simple, sensitive, robust assay for the quantification of carbohydrates even when present in low concentration. The assay can be used for the quantitative determination of any carbohydrates that contain neutral sugars and uronic acids. The reaction is a modification of phenol-sulfuric acid method.</p> <p>In this assay, carbohydrates are hydrolyzed and converted to hydroxyfurfural and furfural, which then react with a highly sensitive developer to form a colored complex, detected spectrophotometrically at 500 nm.</p>
<b>Notes</b>	<p>This product is manufactured by BioVision, an Abcam company and was previously called K2049 High Sensitivity Carbohydrate Assay Kit (Colorimetric). K2049-100 is the same size as the 100 test size of ab284543.</p>
<b>Platform</b>	Microplate (12 x 8 well strips)

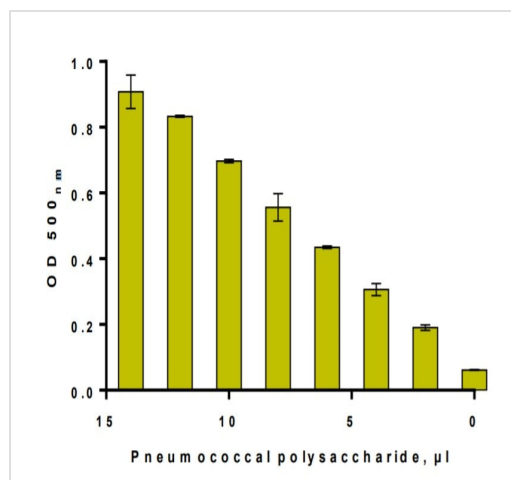
### Properties

---

**Storage instructions** Store at +4°C. Please refer to protocols.

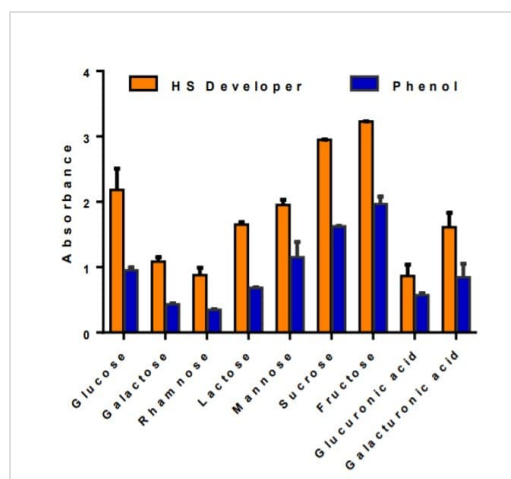
Components	100 tests
Carbohydrate Developer	1 x 1.5ml
Standard (D-Glucose, 2 mg/ml)	1 x 0.2ml

## Images



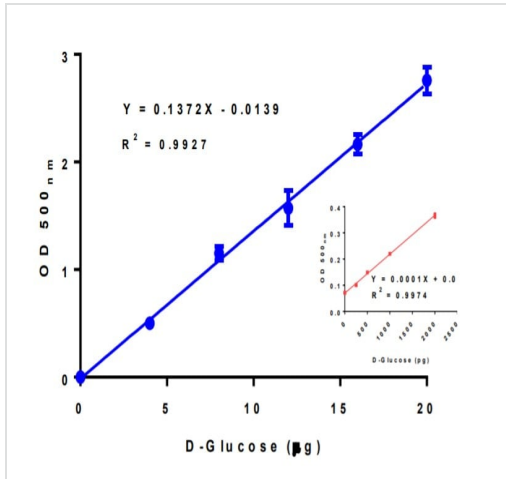
Total Polysaccharide content in Pneumococcal Polysaccharide Powder Type 4 (US Type 4)

Total Polysaccharide content in Pneumococcal Polysaccharide Powder Type 4 (US Type 4), calculated total carbohydrate concentration (glucose equivalent) is  $0.47 \pm 0.02 \mu\text{g}/\mu\text{L}$ . The assay was performed according to the kit assay protocol.



Absorbance of Monosaccharides

Absorbance of monosaccharides upon reacting with High Sensitivity (HS) Carbohydrate Developer and Phenol. Assays were performed following the kit protocol.



D-Glucose Standard Curve. Assays were performed following the kit protocol.

D-Glucose Standard Curve

**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