

Total Carbohydrate Assay Kit - Quantification ab155891

★★★★★ [2 Abreviews](#) [6 References](#) [画像数 4](#)

製品の概要

製品名	Total Carbohydrate Assay Kit - Quantification
検出方法	Colorimetric
サンプルの種類	Tissue, Adherent cells, Suspension cells, Tissue Extracts
アッセイタイプ	Quantitative
全工程の試験時間	0h 30m
製品の概要	Total Carbohydrate Assay (ab155891) is a simple, sensitive and robust method of detecting virtually all carbohydrates.

The carbohydrate assay protocol is based on the phenol - sulfuric acid assay. In the assay, polysaccharides (mono, di, tri, etc.) and their derivatives, in the presence of sulfuric acid, are hydrolyzed to monomers and converted to furfural or hydroxyfurfural. These react with the developer to form a chromogen that can be quantified by measuring the absorbance at $\lambda = 490$ nm.

The total carbohydrate assay kit can detect most forms of carbohydrates, including simple and complex saccharides, glycans, glycoproteins and glycolipids.

Total carbohydrate assay protocol summary:

- add samples and glucose standards to wells
- add concentrated sulfuric acid and incubate for 15 min at 90°C
- add developer and incubate for 5 min
- analyze with a microplate reader

特記事項	<p>This product is manufactured by BioVision, an Abcam company and was previously called K645 Total Carbohydrate Colorimetric Assay Kit. K645-100 is the same size as the 100 test size of ab155891.</p> <p>Carbohydrates play important structural as well as chemical roles in all living systems. Detection of total carbohydrates, therefore, has wide applications.</p>
------	--

試験プラットフォーム	Microplate reader
------------	-------------------

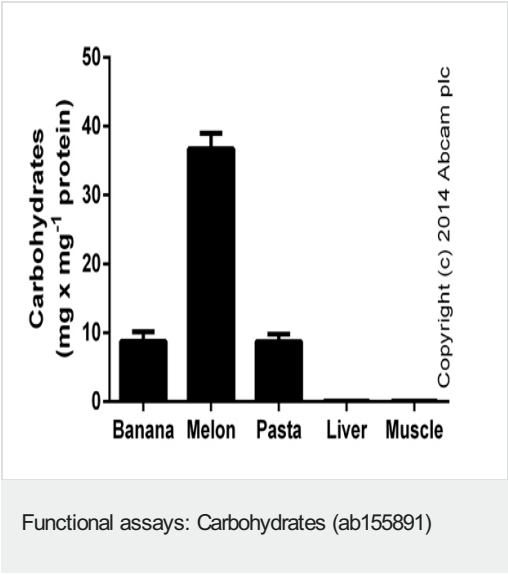
製品の特性

保存方法

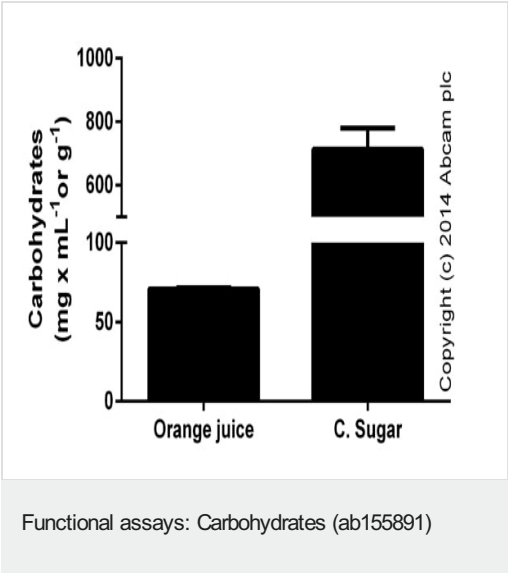
Store at +4°C. Please refer to protocols.

内容	100 tests
Assay Buffer VII	1 x 25ml
Developer Solution IV	1 x 3ml
D-Glucose Standard	1 x 0.2ml

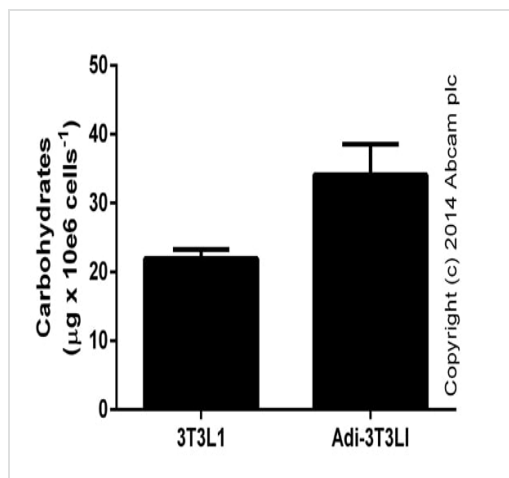
画像



Carbohydrates in food items and mouse tissue lysates, shown as mg per mg of extracted protein (duplicates; +/- SD).

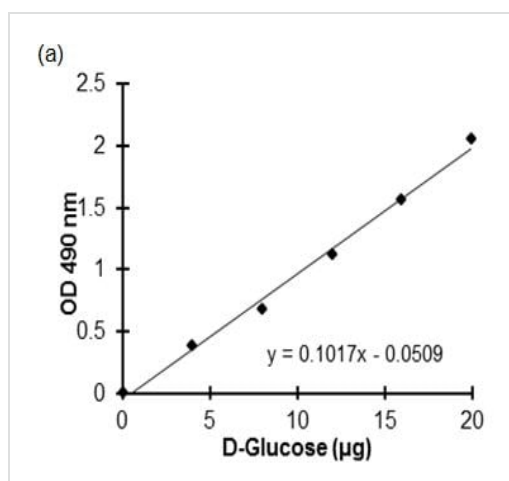


Carbohydrates measured in orange juice and caster sugar. Background signal subtracted, duplicates; +/- SD.



Functional assays: Carbohydrates (ab155891)

Carbohydrates from lysed non-differentiated or adipose differentiated 3T3L1 cells (duplicates; +/- SD). Cells were differentiated using a mixture of small molecules: [ab120840](#), [ab120743](#) and [ab123768](#).



D-Glucose standard curve

D-Glucose standard curve generated using the method described in the protocol booklet.

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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