

Biotin Assay Kit (Colorimetric) ab185441

[2 References](#) [2 Images](#)

Overview

Product name	Biotin Assay Kit (Colorimetric)
Detection method	Colorimetric
Sample type	Purified protein
Assay type	Semi-quantitative
Sensitivity	15 ng
Product overview	Abcam's Biotin Quantitation Kit (Colorimetric) (ab185441) is based on the differential binding of Streptavidin to Biotin and a dye, 2-(4-hydroxyazobenzene) benzoic acid (HABA). When HABA is bound to Streptavidin, addition of free biotin or biotinylated biomolecules results in its displacement in the reaction. These changes are manifested in terms of decrease in the overall absorption (OD 500 nm) and can be quantitatively correlated to the amount of biotin present in solution. The kit also provides Biotin as a Standard and Biotinylated BSA as a positive control. This kit can be used to estimate as low as 60 pmol or 15 ng of biotin in solution.

Notes

This product is manufactured by BioVision, an Abcam company and was previously called K811 Biotin Quantitation Kit (Colorimetric). K811-100 is the same size as the 100 test size of ab185441.

Biotin (Vitamin H) is an important biomolecule that has been widely used in Biotinylation reactions to label antibodies or other proteins of interest. Due to its low molecular weight, biotin usually does not cause any significant changes in protein conformation and biological activity. Biotinylated proteins/antibodies are widely isolated and assayed with streptavidin pull-down assay, affinity chromatography, ELISA and Western blotting etc. It is often desired to determine the degree of substitution of such biotinylated biomolecules.

Platform Microplate reader

Properties

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

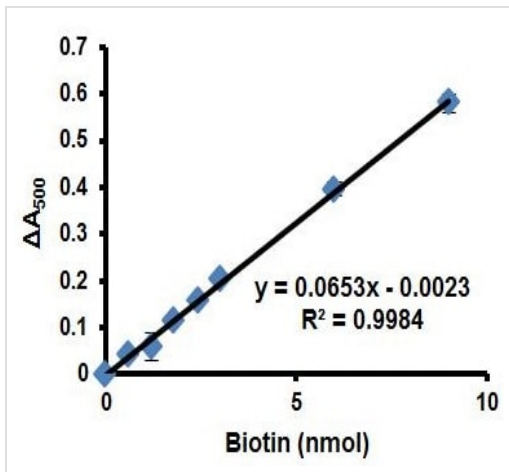
Components	100 tests
Biotin Assay Buffer	1 x 25ml
Biotin Assay Reagent A	1 x 8ml

Components	100 tests
Biotin Assay Reagent B	1 x 1ml
Biotin Standard	1 x 100µl
Biotinylated BSA	1 x 200µl

Relevance

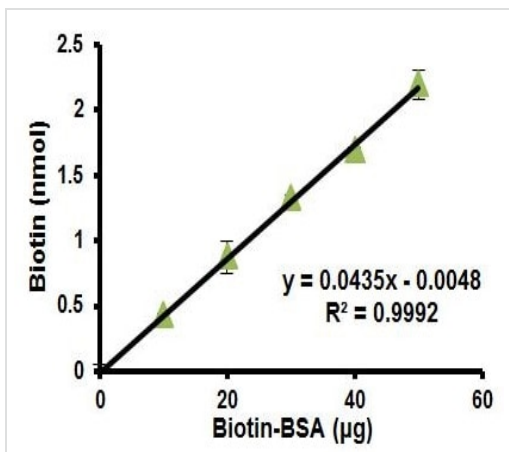
Biotin is a water soluble vitamin, generally classified as a B complex vitamin, also called vitamin B4. After the initial discovery of biotin, nearly forty years of research were required to establish it as a vitamin. Biotin is required by all organisms but can only be synthesized by bacteria, yeasts, molds, algae, and some plant species. Biotin is required as prosthetic group of enzymes involved in incorporation of carbon dioxide into organic compounds. Biotin has a MW of 244 Da.

Images



This is example data only.

Biotin Standard Curve



Linear plot obtained for the amount of biotin (nmol) vs amount of biotinylated BSA. Assays were performed according to the kit protocol. This is example data only.

Amount of biotin vs amount of biotinylated BSA

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