

Bradford Assay Kit ab102535

25 References **画像数 3**

製品の概要

製品名	Bradford Assay Kit
検出方法	Colorimetric
サンプルの種類	Serum, Plasma, Cell culture extracts, Other biological fluids, Tissue Extracts
アッセイタイプ	Quantitative
全工程の試験時間	0h 20m
製品の概要	<p>Bradford Assay Kit ab102535 provides a simple and rapid procedure for determining the concentration of protein in solution.</p> <p>The Bradford assay protocol uses an improved Coomassie blue G reagent which forms a blue complex in the presence of protein. The intensity of the blue complex is proportional to the amount of protein in the sample and can be easily measured by spectrophotometer or plate reader at 595 nm.</p> <p>Bradford assay protocol summary:</p> <ul style="list-style-type: none"> - prepare serial dilutions of protein standard - add samples and standards to wells or tubes - add Bradford protein assay solution and incubate for 5 min at room temp - measure absorbance at 595 nm <p>Support on how to determine protein concentration can be found here</p>
特記事項	This product is manufactured by BioVision, an Abcam company and was previously called K810 Protein Quantitation Kit. K810-1000 is the same size as the 1000 test size of ab102535.
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.

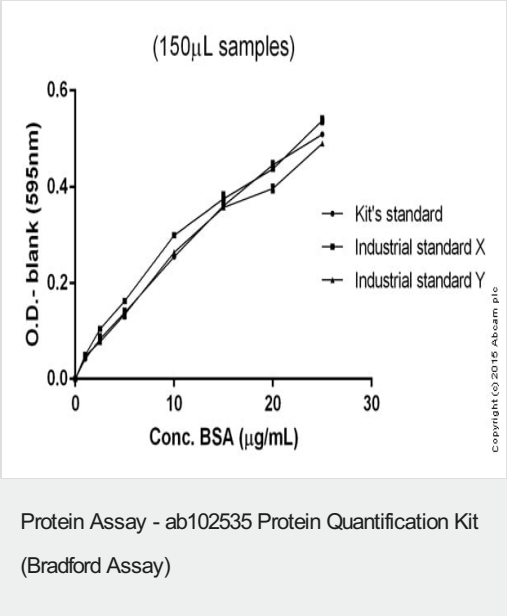
内容	1000 tests
5X Protein Assay Solution	1 x 20ml

内容	1000 tests
Protein Standard I	1 x 1.5ml

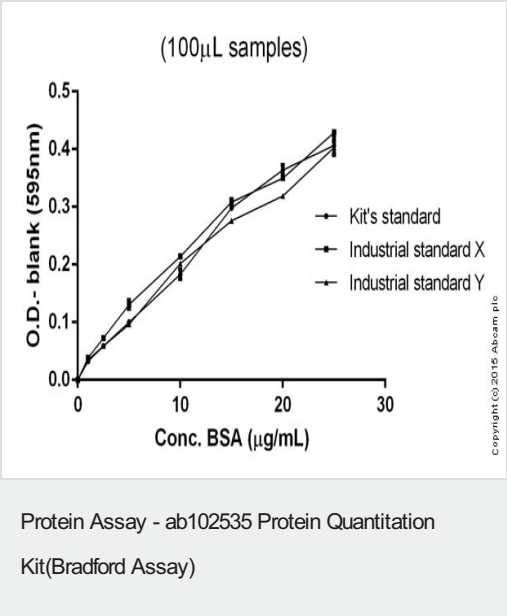
関連性

Protein Quantitation is a procedure for determining the concentration of protein in solution.

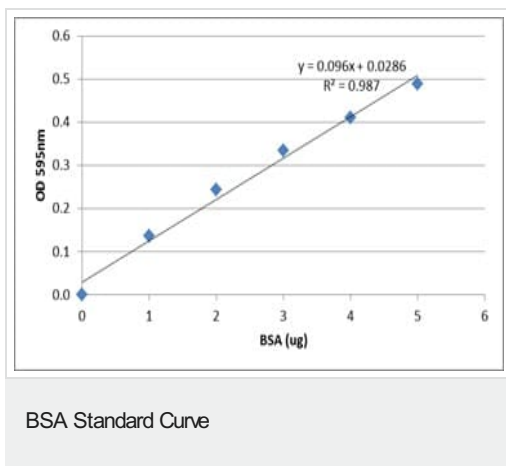
画像



Colorimetric standard curves of 150 microliter samples: mean of duplicates with background readings subtracted.



Colorimetric standard curves of 100 microliter samples: mean of duplicates with background readings subtracted.



Colorimetric standard curve of 10 microliter samples: mean of duplicates with background readings subtracted.

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