

Product datasheet

Protein Quantitation Kit (Bradford Assay) ab102535

3 References [画像数 3](#)

製品の概要

製品名	Protein Quantitation Kit (Bradford Assay)
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	Abcam's Protein Quantitation Kit (Bradford Assay) provides a simple and rapid procedure for determining the concentration of protein in solution. The method utilizes 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.
アプリケーション	適用あり: Functional Studies

法規制情報

医薬用外毒物

製品の特性

保存方法 Store at +4°C. Please refer to protocols.

内容	ラベル	1000 tests
Protein Assay Solution	NM	1 x 20ml
Protein Standard	Yellow	1 x 1.5ml

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

アプリケーション

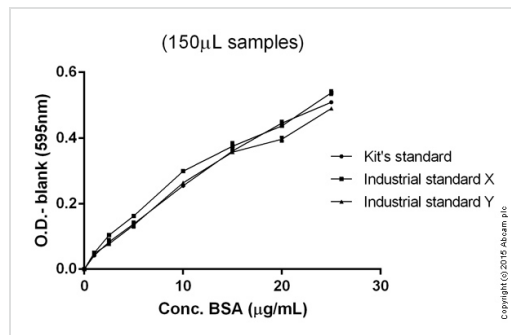
Our [Abpromise guarantee](#) covers the use of **ab102535** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Functional Studies

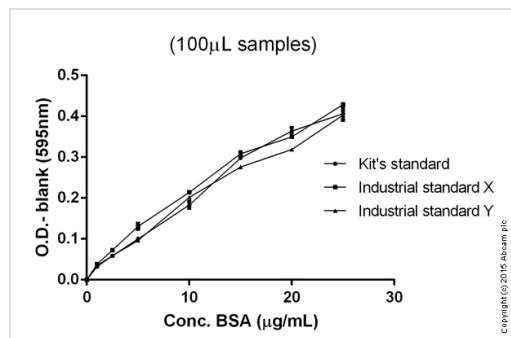
Use at an assay dependent dilution.

画像



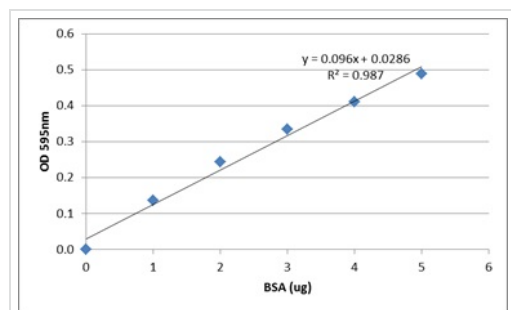
Colorimetric standard curves of 150 microliter samples: mean of duplicates with background readings subtracted.

Protein Assay - ab102535 Protein Quantification Kit (Bradford Assay)



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

Protein Assay - ab102535 Protein Quantitation Kit(Bradford Assay)



BSA Standard Curve

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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 <http://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