

Griess reagent ab234044

5 References [画像数 2](#)

医薬用外劇物

製品の概要

製品名	Griess reagent
検出方法	Colorimetric
サンプルの種類	Urine, Serum, Plasma, Cell Lysate, Tissue Lysate
アッセイタイプ	Quantitative
検出感度	1 nmol/well
全工程の試験時間	1h 30m
製品の概要	Griess Reagent Kit ab234044 uses a classic protocol for the estimation of nitrite in biological samples.

Nitrite is reduced to nitrogen oxide using Griess Reagent I. Nitrogen oxide then reacts with Griess Reagent II forming a stable product that can be detected by its absorbance at 540 nm.

The two-step assay is simple, fast and can detect nitrite levels as low as 1 nmol/well.

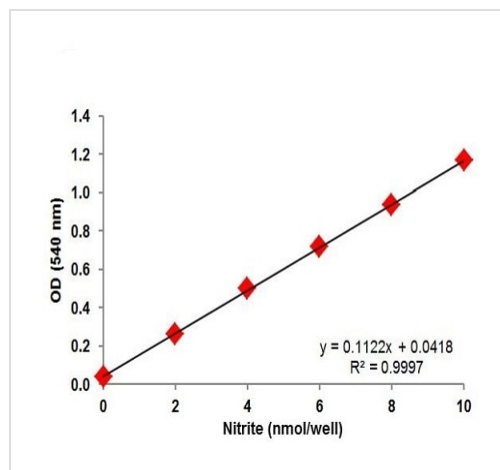
特記事項 This product is manufactured by BioVision, an Abcam company and was previously called K544 Nitrite Assay Kit (Griess Reagent). K544-200 is the same size as the 200 test size of ab234044.

試験プラットフォーム Microplate reader

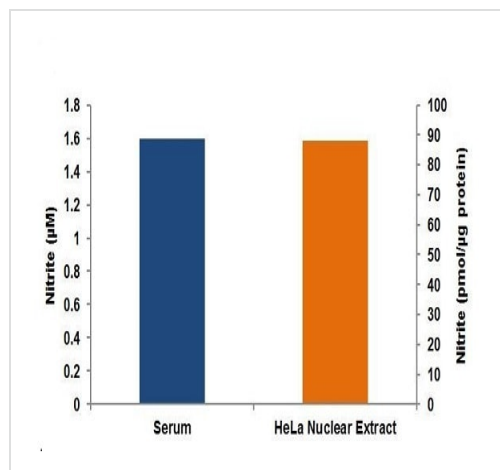
製品の特性

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

内容	200 tests	500 tests
Buffer III	1 x 30ml	1 x 75ml
Griess Reagent I	1 x 10ml	1 x 25ml
Griess Reagent II	1 x 10ml	1 x 25ml
Nitrite Standard	1 vial	1 vial



Nitrite standard curve



Nitrite concentration in pooled human serum and HeLa nuclear extracts

Human serum was deproteinized using 10 kDa columns. Filtrate (50 µl, undiluted) was collected and assayed according to kit protocols. HeLa nuclear cells were homogenized using Nitrite Assay Buffer (100 µg protein).

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