

Nitric Oxide Assay Kit ab272517

5 References [画像数 1](#)

医薬用外劇物

製品の概要

製品名	Nitric Oxide Assay Kit
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Saliva, Urine, Serum, Plasma, Whole Blood, Tissue Extracts, Cell Lysate
アッセイタイプ	Quantitative
検出範囲	0.6 μ M - 200 μ M
製品の概要	<p>Nitric Oxide Assay Kit (ab272517) is a simple, direct and automation-ready procedure for measuring NO. Since NO is oxidized to nitrite and nitrate, it is common practice to quantitate total $\text{NO}_2^-/\text{NO}_3^-$ as a measure for NO level. This Nitric Oxide Assay Kit is designed to accurately measure NO production following reduction of nitrate to nitrite using an improved Griess method. The procedure is simple and the time required for sample pre-treatment and assay is reduced to as little as 30 minutes.</p> <p>Sensitive and accurate: Detection range 0.6 - 200 μM in 96-well plate.</p> <p>Rapid and reliable: Using an optimized VCl_3 reagent, the time required for reduction of NO_3^- to NO_2^- is 10 min at 60°C.</p> <p>Simple and high-throughput: The procedure involves mixing sample with three reagents, incubation for 10 minutes at 60°C and reading the optical density. Can be readily automated to measure thousands of samples per day.</p>
アプリケーション	適用あり: Functional Studies
試験プラットフォーム	Microplate

製品の特性

保存方法 Store at -20°C. Please refer to protocols.

内容	100 tests
NaOH	1 x 1ml

内容	100 tests
Nitrite Standard (1 mM)	1 x 1ml
Reagent A	1 x 12ml
Reagent B	1 x 500µl
Reagent C	1 x 12ml
ZnSO4	1 x 1ml

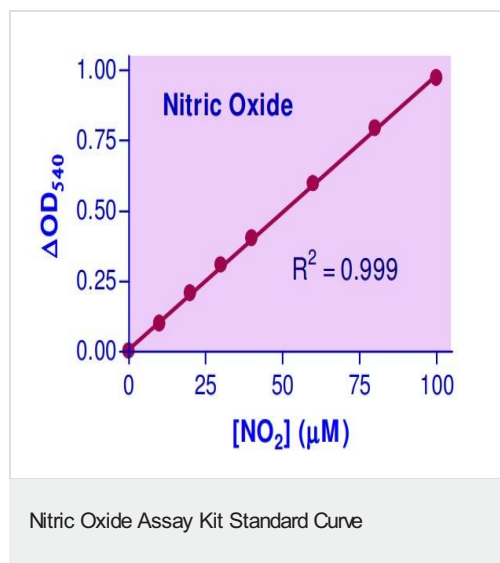
関連性 Nitric oxide (NO) is a key vertebrate biological messenger, playing an important role in neurotransmission, vascular regulation, immune responses and apoptosis. NO, also known as "endothelium-derived relaxing factor" or "EDRF", is synthesized from L-arginine, oxygen and NADPH by various NO synthases. Most of the NO in the cell is oxidized to nitrite (NO_2^-) and nitrate (NO_3^-), and therefore the concentrations of these anions are generally as a quantitative measure of NO production.

アプリケーション

The Abpromise guarantee Abpromise保証は、 次のテスト済みアプリケーションにおけるab272517の使用に適用されます
 アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご検討ください。

アプリケーション	Abreviews	特記事項
Functional Studies		Use at an assay dependent concentration.

画像



Typical standard curve – data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.

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