

Ammonia Assay Kit ab272538

画像数 1

医薬用外劇物

製品の概要

製品名	Ammonia Assay Kit
検出方法	Fluorescent
サンプルの種類	Urine, Other biological fluids
アッセイタイプ	Quantitative
検出範囲	0.012 mM - 1 mM
製品の概要	<p>Ammonia Assay Kit (ab272538) is simple, direct and automation-ready procedures for measuring NH₃ are very desirable. This assay is based on an improved o-phthalaldehyde method. This reagent reacts with ammonia/ ammonium and forms a fluorescent product. The fluorescence intensity ($\lambda_{\text{ex/em}}$ = 360/450nm) is proportional to the ammonia concentration in the sample.</p> <p>Fast and sensitive: Linear detection range of 0.012 - 1 mM ammonia.</p> <p>Convenient and high-throughput: Homogeneous "mix-incubate-measure" type assay. Can be readily automated on HTS liquid handling systems for processing 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	200 tests
Assay Buffer	1 x 20ml
Reagent A	1 x 1ml
Reagent B	1 x 1ml
Standard (20mM Ammonium chloride)	1 x 400µl

関連性

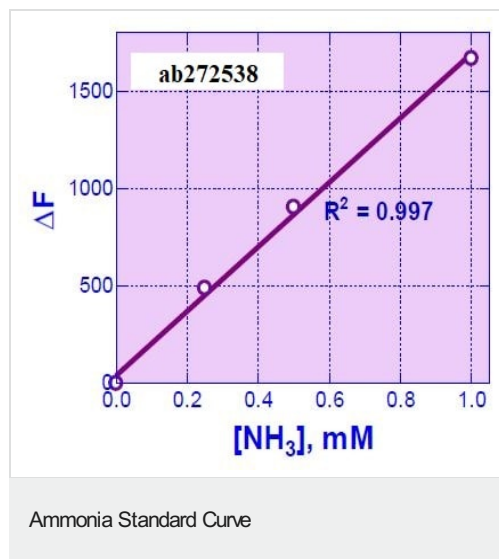
Ammonia is an important source of nitrogen for living systems. Nitrogen is required for the synthesis of amino acids, which are the building blocks of protein. Ammonia is a metabolic product which is created through amino acid deamination. It plays an important role in both normal and abnormal animal physiology like normal animal acid/base balance.

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ab272538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確認ください。

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