

Product datasheet

Ammonia Assay Kit ab272538

1 Image

Overview

| | |
|----------------------------|---|
| Product name | Ammonia Assay Kit |
| Detection method | Fluorescent |
| Sample type | Urine, Other biological fluids |
| Assay type | Quantitative |
| Range | 0.012 mM - 1 mM |
| Product overview | <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 ($\lambda_{ex/em} = 360/450\text{nm}$) 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> |
| Tested applications | Suitable for: Functional Studies |
| Platform | Microplate |

Properties

Storage instructions Store at -20°C. Please refer to protocols.

| Components | 200 tests |
|-----------------------------------|-----------|
| Assay Buffer | 1 x 20ml |
| Reagent A | 1 x 1ml |
| Reagent B | 1 x 1ml |
| Standard (20mM Ammonium chloride) | 1 x 400µl |

Relevance

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.

Applications

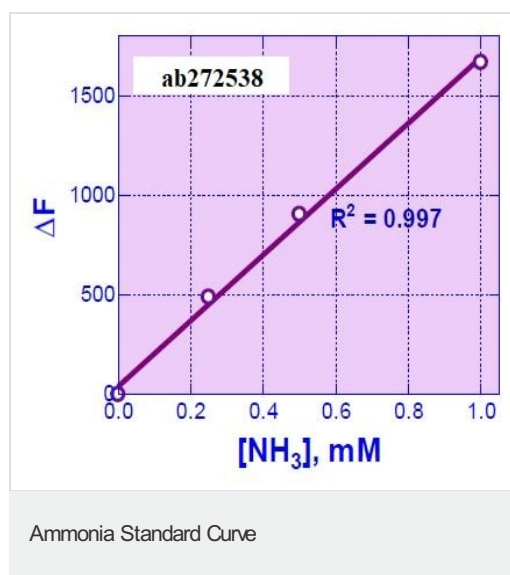
The Abpromise guarantee

Our [Abpromise guarantee](#) covers the use of ab272538 in the following tested applications.

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

| Application | Abreviews | Notes |
|--------------------|-----------|--|
| Functional Studies | | Use at an assay dependent concentration. |

Images



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.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors