

Product datasheet

Ammonia Assay Kit - Modified Berthelot - (Colorimetric) ab102509

[5 References](#) [2 Images](#)

Overview

Product name	Ammonia Assay Kit - Modified Berthelot - (Colorimetric)
Sample type	Urine, Serum, Plasma, Other biological fluids, Cell culture media
Assay type	Quantitative
Sensitivity	< 10 μ M
Assay time	0h 30m
Product overview	<p>Ammonia Assay Kit - Modified Berthelot (Colorimetric) ab102509 provides a rapid, simple, sensitive, and reliable assay to measure ammonia.</p> <p>In this non-enzymatic ammonia assay protocol, ammonia is used to form indophenol, a highly colored product easily quantifiable by colorimetry ($\lambda=670\text{nm}$) using a plate reader.</p> <p>The kit can detect less than 1 nmol($\sim 10 \mu\text{M}$) ammonia or ammonium using either a 96 well or 384 well microwell plate.</p> <p>Ammonia assay protocol summary:</p> <ul style="list-style-type: none"> - add samples and standards to wells - add reagent 1 and reagent 2 and incubate for 30 min - analyze with a microplate reader
Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K470 Ammonia Colorimetric Assay Kit II. K470-100 is the same size as the 100 test size of ab102509.</p> <p><u>Ammonia assay kit ab83360</u> is an alternative to this product, and is the most popular ammonia assay kit.</p>
Platform	Microplate reader

Properties

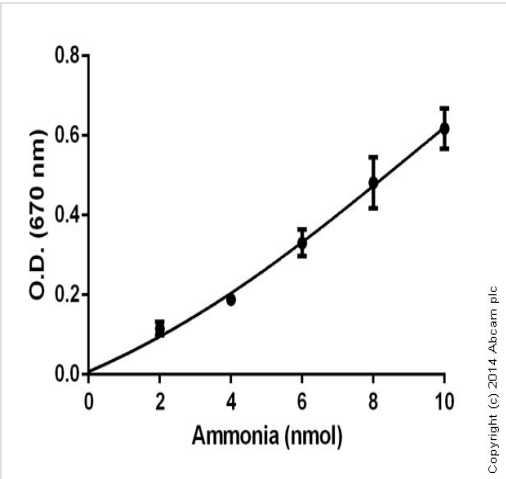
Storage instructions	Store at +4°C. Please refer to protocols.
-----------------------------	---

Components	100 tests
Reagent II	1 x 8ml
Reagent III	1 x 4ml
Ammonium Standard II	1 x 0.1ml

Relevance

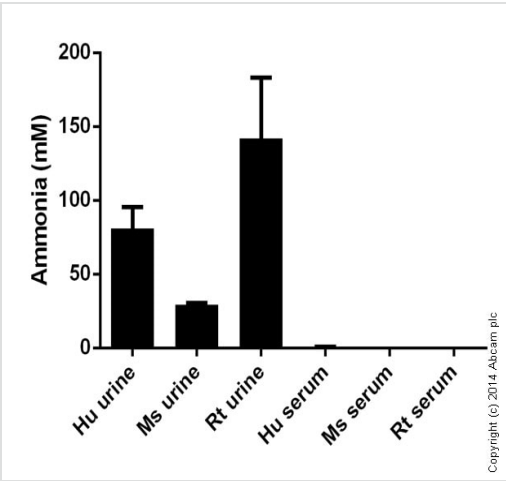
Ammonia is an important source of nitrogen for living systems. Nitrogen is required for the synthesis of amino acids, which are the building blocks of protein. Ammonia is a metabolic product which is created through amino acid deamination. It plays an important role in both normal and abnormal animal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

Images



Standard curve: mean of duplicates (+/- SD) with background reads subtracted.

Functional Studies - Ammonia Colorimetric Assay
Kit (Modified Berthelot Assay) (ab102509)



Ammonia measured in biological fluids showing quantity (mM) in tested samples. Samples were diluted 10-1000 fold.

Functional Studies - Ammonia Colorimetric Assay
Kit (Modified Berthelot Assay) (ab102509)

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