

Product datasheet

Ammonia Assay Kit - Modified Berthelot - (Colorimetric) ab102509

2 图像

概述

产品名称	Ammonia Assay试剂盒- Modified Berthelot - (Colorimetric)
样品类型	Urine, Serum, Plasma, Other biological fluids, Cell culture media
检测类型	Quantitative
灵敏度	< 10 μ M
检测时间	0h 30m
产品概述	Abcam's Ammonia Assay Kit - Modified Berthelot - (Colorimetric) (ab102509) provides a rapid, simple, sensitive, and reliable assay suitable for research and high throughput assay of Ammonia. In this non-enzymatic assay, ammonia is used to form indophenol, a highly colored product easily quantifiable by colorimetry ($\lambda=670\text{nm}$) using a plate reader. Visit our FAQs page for tips and troubleshooting.

说明 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.

Store the kit at +4°C.

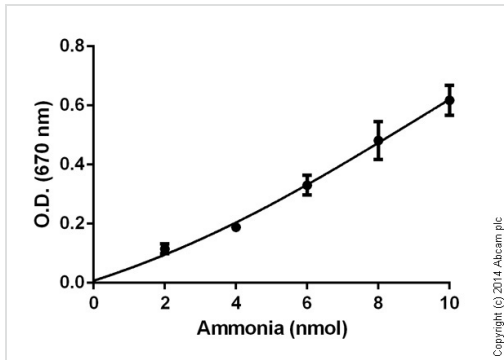
性能

存放说明 Store at +4°C. Please refer to protocols.

组件	标识符	100 tests
Ammonia Reagent 1	Amber	1 x 8ml
Ammonia Reagent 2	Clear	1 x 4ml
Ammonium Chloride Standard	Yellow	1 x 0.1ml

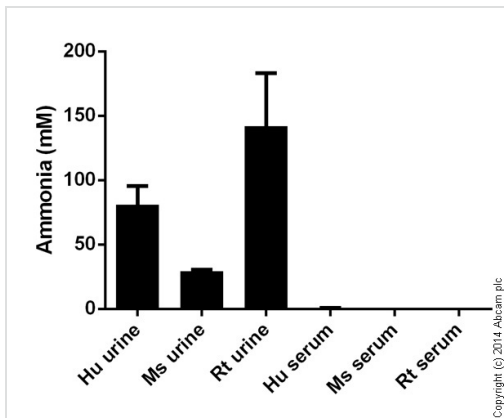
相关性 Ammonia is an important source of nitrogen for living systems. Nitrogen is required for the synthesis of amino acids, which are the building blocks of protein. Ammonia is a metabolic product which is created through amino acid deamination. It plays an important role in both normal and abnormal animal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

图片



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 AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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.cn/abpromise> or contact our technical team.

Terms and conditions

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