

## Product datasheet

# Ammonia Assay Kit - Modified Berthelot - (Colorimetric) ab102509

### 2 图像

#### 概述

<b>产品名称</b>	Ammonia Assay试剂盒- Modified Berthelot - (Colorimetric)
<b>样品类型</b>	Urine, Serum, Plasma, Other biological fluids, Cell culture media
<b>检测类型</b>	Quantitative
<b>灵敏度</b>	< 10 $\mu$ M
<b>检测时间</b>	0h 30m
<b>产品概述</b>	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 <a href="#">FAQs page</a> for tips and troubleshooting.
<b>说明</b>	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.
<b>经测试应用</b>	<b>适用于:</b> Functional Studies

#### 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.

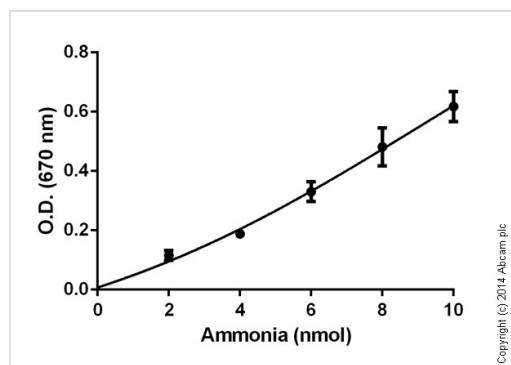
## 应用

Our [Abpromise guarantee](#) covers the use of **ab102509** in the following tested applications.

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

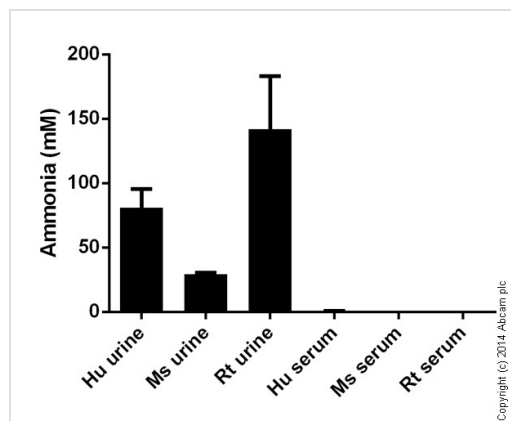
应用	Ab评论	说明
Functional Studies		Use at an assay dependent dilution.

## 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 <http://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