

Product datasheet

Formaldehyde Assay Kit ab272524

1 Image

Overview

Product name	Formaldehyde Assay Kit
Detection method	Fluorescent
Sample type	Urine, Serum, Plasma, Other biological fluids, Cell Lysate
Assay type	Quantitative
Sensitivity	1.5 µM
Product overview	<p>Formaldehyde Assay Kit (ab272524) provides a convenient fluorimetric means to measure formaldehyde in biological samples. In the assay, formaldehyde is derivatized with acetoacetanilide in the presence of ammonia. The resulting fluorescent product is then quantified fluorimetrically ($\lambda_{exc}/\lambda_{em} = 370/470nm$). The assay is simple, sensitive, stable and high-throughput adaptable. The assay can detect as low as 1.5 µM formaldehyde in biological samples.</p> <p>Safe. Non-radioactive assay.</p> <p>Sensitive and accurate. As low as 1.5 µM (45 ppb) formaldehyde can be quantified.</p> <p>Homogeneous and convenient. "Mix-incubate-measure" type assay. No wash and reagent transfer steps are involved.</p> <p>Robust and amenable to HTS: 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 +4°C. Please refer to protocols.

Components	100 tests
Neutralizer	2 x 1.5ml

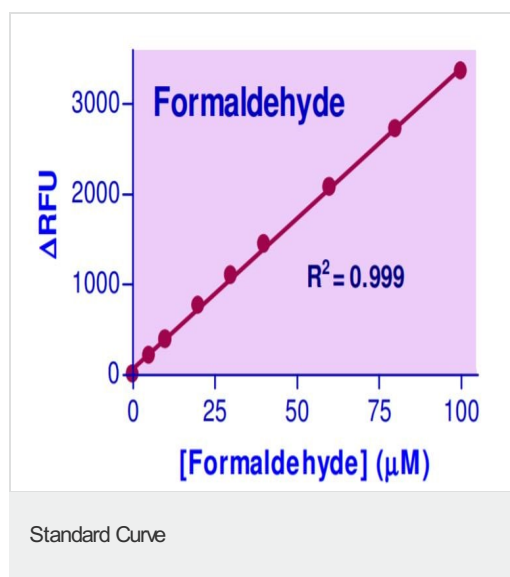
Components	100 tests
Reagent A	1 x 5ml
Reagent B	1 x 3ml
Standard	1 x 100µl
TCA (10%)	1 x 5ml

Applications

The Abpromise guarantee Our [Abpromise guarantee](#) covers the use of ab272524 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