

Product datasheet

Free Amino Nitrogen Assay Kit ab272523

1 Image

Overview

Product name	Free Amino Nitrogen Assay Kit
Detection method	Colorimetric
Sample type	Food samples
Assay type	Quantitative
Range	0.2 mM - 10 mM
Product overview	<p>Free Amino Nitrogen Assay Kit (ab272523) measures alpha amino acids, ammonia, and end group amino nitrogens. The ninhydrin based reaction is a superior method for determining only alpha amino acids and ammonia compared to the traditional Kjeldahl, which measures nitrogen from all sources. Only requiring low sample volumes, the stable ninhydrin reagent provides a simple and accurate method for determining Free Amino Nitrogen concentrations.</p> <p>Fast and sensitive. Linear detection range (5 µL sample): 0.2 to 10 mM for 10 min reaction.</p> <p>Convenient and high-throughput. Homogeneous "mix-incubate-measure" type assay. Can be readily automated to process 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
Glycine Standard (20 mM)	1 x 500µl
Reagent A	1 x 18ml
Reagent B	1 x 600µl

Applications

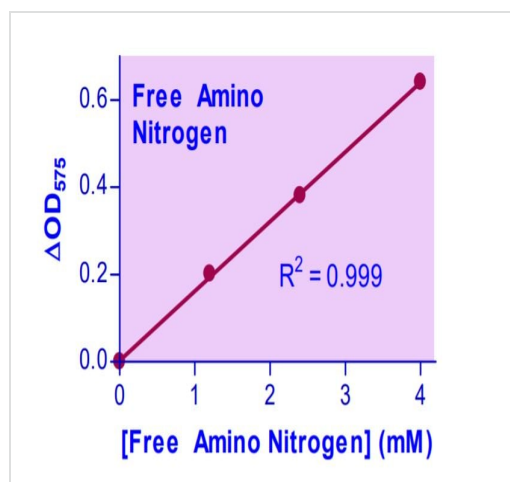
The Abpromise guarantee

Our [Abpromise guarantee](#) covers the use of ab272523 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



Free Amino Nitrogen Assay Kit Standard Curve

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