

# L-Amino Acid Assay Kit ab65347

**22 References**    **画像数 3**

### 製品の概要

製品名	L-Amino Acid Assay Kit
検出方法	Colorimetric/Fluorometric
サンプルの種類	Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids, Tissue Extracts
アッセイタイプ	Quantitative
検出感度	$\leq 40 \mu\text{M}$
全工程の試験時間	0h 30m
製品の概要	L-Amino Acid Assay Kit ab65347 is used measure the level of L-amino acids in biological samples. There is no requirement for sample pretreatment or purification.

The L-amino acid(s) level can be quantified using fluorometric (at Ex/Em = 535/587 nm) or colorimetric (at  $\lambda = 570 \text{ nm}$ ) methods in 96-well plates.

Amino acid assay protocol summary:

- add reaction mix to sample and standard wells
- incubate for 30 min
- analyze with microplate reader

特記事項	<p>This product is manufactured by BioVision, an Abcam company and was previously called K639 L-Amino Acid Quantitation Colorimetric/Fluorometric Kit. K639-100 is the same size as the 100 test size of ab65347.</p> <p>L-Amino acids are the most essential elements in biology. Accurately quantitating L-amino acids in body fluids or purified samples may provide valuable information for diagnostic or basic research studies.</p>
------	--

試験プラットフォーム	Microplate reader
------------	-------------------

### 製品の特性

保存方法	Store at -20°C. Please refer to protocols.
------	--

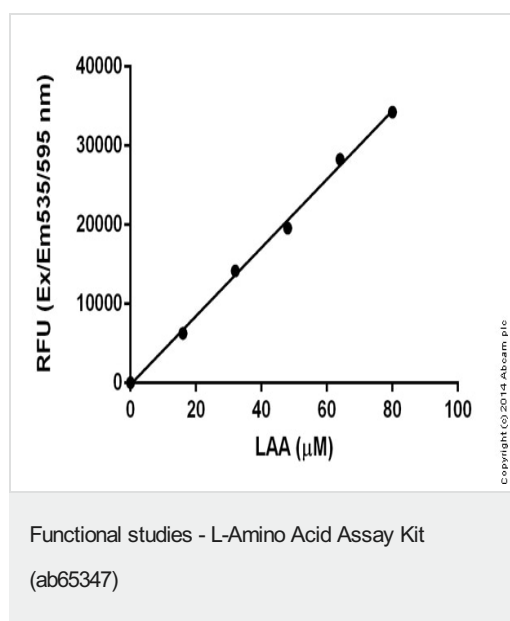
内容	100 tests
Assay Buffer II	1 x 25ml

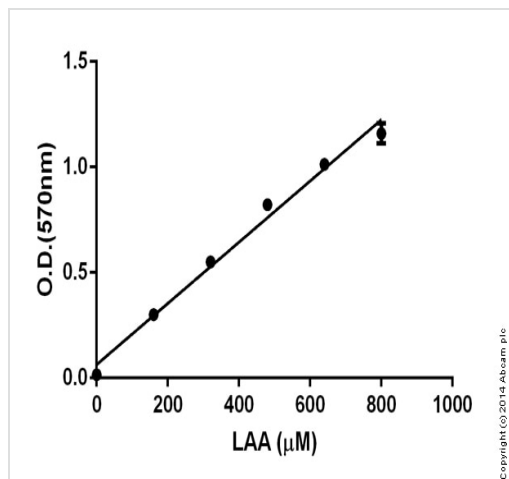
内容	100 tests
L-Amino Acid Enzyme Mix	1 vial
L-Amino Acid Standard	1 x 300μl
OxiRed Probe	1 x 0.2ml

**関連性**

Amino acids perform various important functions for the body and serve as the materials for the body's cells, hormones, and enzymes. Of the standard amino acids, all but glycine can exist in either of the two optical isomers, called D and L enantiomers, which are mirror images of each other. The L-amino acids represent the vast majority of amino acids found in proteins during translation in the ribosome, and are therefore the most essential element in biology and metabolic functions.

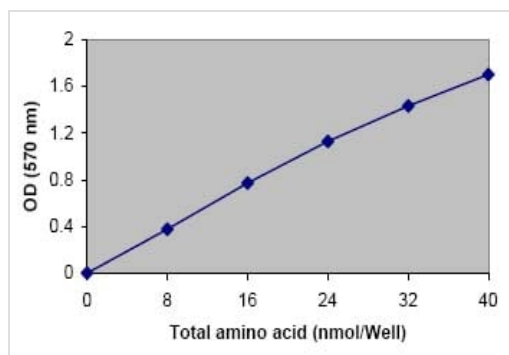
## 画像





Standard curve (colorimetric): mean of duplicates (+/-SD) with background readings subtracted

Functional studies - L-Amino Acid Assay Kit  
(ab65347)



L-Amino Acid Standard Curve.

Functional Studies - L-Amino Acid Assay Kit  
(ab65347)

**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.co.jp/abpromise> or contact our technical team.

### Terms and conditions

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