

Asparaginase Activity Assay Kit ab107922

[5 References](#) [4 Images](#)

Overview

Product name	Asparaginase Activity Assay Kit
Detection method	Colorimetric/Fluorometric
Sample type	Cell Lysate
Assay type	Enzyme activity
Assay time	0h 40m
Product overview	Asparaginase Activity Assay Kit (Colorimetric/Fluorometric) (ab107922) provides a simple, direct and automation-ready procedure for measuring asparaginase activity in biological samples.

In the asparaginase assay protocol, asparaginase hydrolyzes asparagine to generate aspartic acid, which can be detected indirectly colorimetrically (OD=570 nm) or fluorescently (Ex/Em = 535/590 nm) using a coupled enzymatic reaction.

Asparaginase assay protocol summary:

- add reaction mix to sample and standard wells, mix and incubate for 10 min
- analyze with microplate reader in kinetic mode for 30-60 min

Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K754 Asparaginase Activity Colorimetric/Fluorometric Assay Kit. K754-100 is the same size as the 100 test size of ab107922.</p> <p>Asparaginase (EC 3.5.1.1) is a homotetramer that catalyzes the hydrolysis of asparagine to aspartic acid and ammonia and exhibits about a 2-4% activity on glutamine and 5% on D-asparagine. Asparaginase does not occur naturally in humans but is found in bacteria, plants and many animals (e.g. guinea pigs).</p>
-------	--

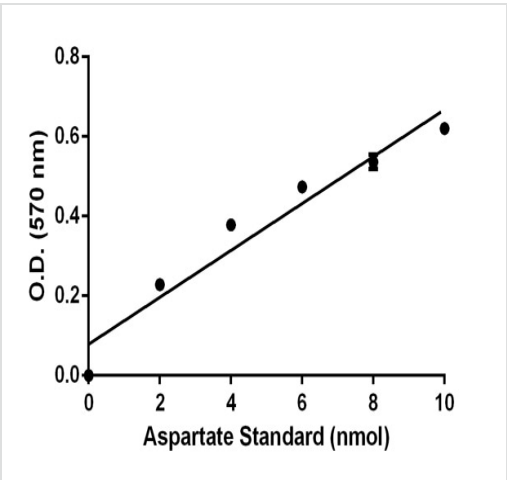
Properties

Storage instructions	Store at -20°C. Please refer to protocols.
----------------------	--

Components	100 tests
Asparaginase Positive Control	1 vial
Aspartate Enzyme Mix	1 vial

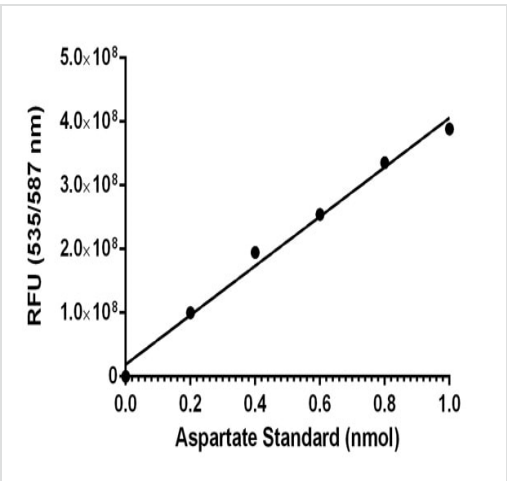
Components	100 tests
Aspartate Standard	1 x 0.1ml
Assay Buffer IV	1 x 25ml
Converter Enzyme XII	1 vial
OxiRed Probe	1 x 0.2ml
Asparaginase Substrate Mix	1 vial

Images



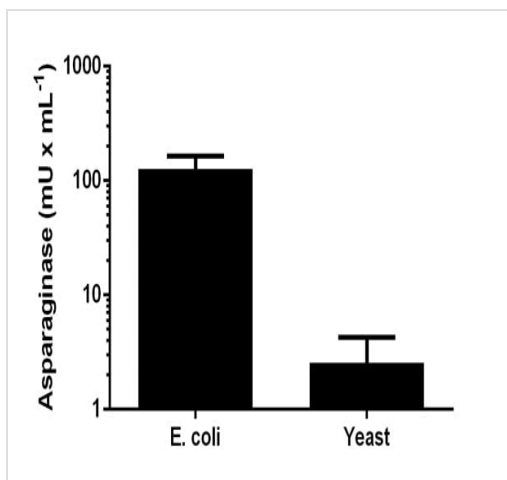
Typical asparaginase standard calibration curve using colorimetric reading.

Functional Studies - Asparaginase Activity Assay Kit (Colorimetric)



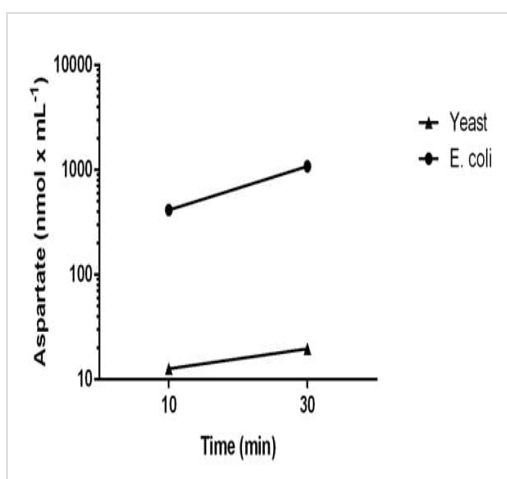
Typical asparaginase standard calibration curve using fluorometric reading.

Functional Studies - Asparaginase Activity Assay Kit (Fluorometric)



Asparaginase measured in cell lysates showing quantity (mU) per mL of tested sample.

Functional Studies - Asparaginase Activity Assay
Kit (Fluorometric)



Aspartate measured in cell lysates after 10 min and 30 min incubation time showing quantity (nmol) per mL of tested sample.

Functional Studies - Asparaginase Activity Assay
Kit (Fluorometric)

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