

Aspartate Assay Kit ab102512

★★★★★ [1 Abreviews](#) [14 References](#) [5 Images](#)

Overview

Product name	Aspartate Assay Kit
Detection method	Colorimetric/Fluorometric
Sample type	Cell culture supernatant, Urine, Serum, Other biological fluids, Tissue Extracts
Assay type	Quantitative
Sensitivity	> 2 µM
Range	2 µM - 200 µM
Assay time	0h 40m
Product overview	<p>Aspartate Assay Kit ab102512 provides a simple, convenient assay to measure aspartate in a variety of samples.</p> <p>In the aspartate assay protocol, aspartate is converted to pyruvate which is oxidized with the conversion of a probe into a highly colored (570 nm) and fluorescent (Ex/Em 535/587 nm) species proportional to the amount of aspartate in samples. Aspartate can be quantified in the range between 0.1–10 nmoles/well (2-200 µM).</p>
Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K552 Aspartate Colorimetric/Fluorometric Assay Kit. K552-100 is the same size as the 100 test size of ab102512.</p> <p>L-Aspartic acid (Asp) is one of the 20 proteinogenic amino acids as building block for protein. Although a non-essential amino acid in mammals, it is a precursor to several other amino acids including four essential amino acids (Met, Thr, Ile and Lys).</p>
Platform	Microplate reader

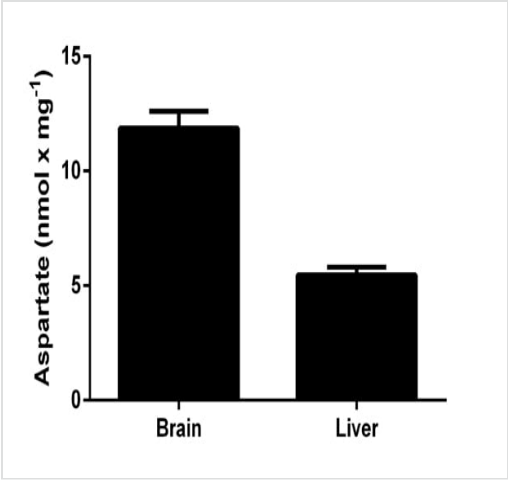
Properties

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

Components	100 tests
Aspartate Enzyme Mix	1 vial
Aspartate Standard	1 x 0.1ml

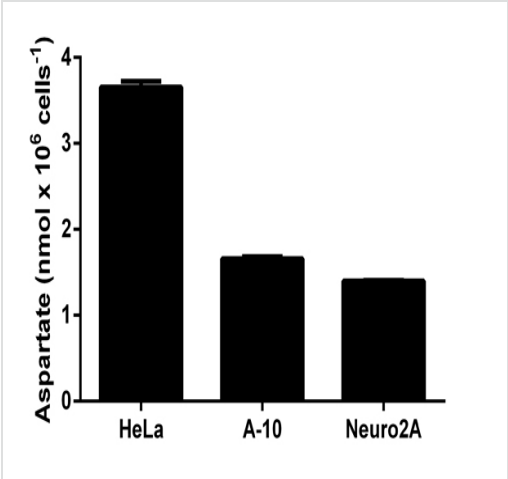
Components	100 tests
Assay Buffer IV	1 x 25ml
Converter Enzyme XII	1 vial
OxiRed Probe	1 x 0.2ml
Sample Clean-Up Mix I	1 vial

Images



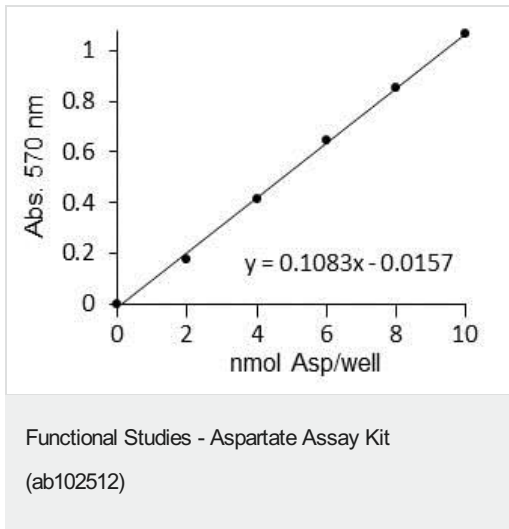
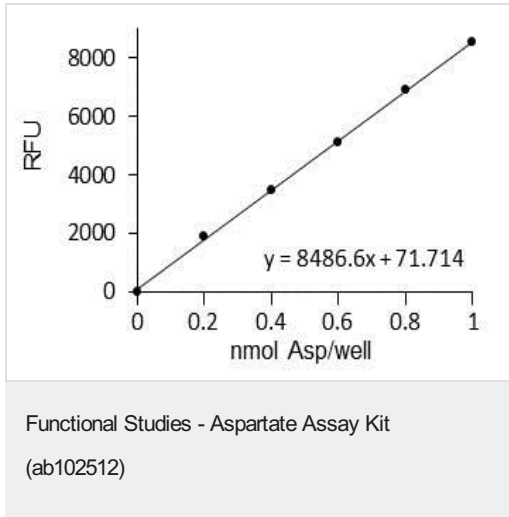
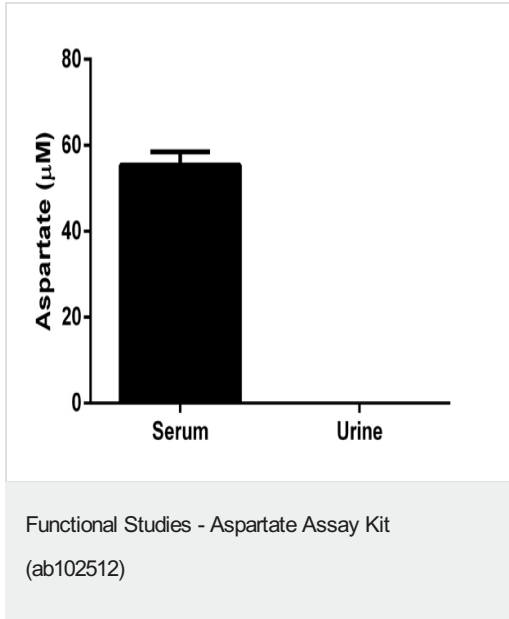
Aspartate levels measured colourimetrically in mouse tissue lysates (mg of extracted protein: background signal subtracted, mean of duplicates; +/- SD).

Functional Studies - Aspartate Assay Kit
(ab102512)



Aspartate levels measured fluorometrically in cell lysates (background signal subtracted, mean of duplicates; +/- SD).

Functional Studies - Aspartate Assay Kit
(ab102512)



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