

Human Aspartate Aminotransferase ELISA Kit ab285299

1 Image

Overview

Product name	Human Aspartate Aminotransferase ELISA Kit			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	Overall			CV%
				< 8%
	Inter-assay			
	Sample	n	Mean	SD
	Overall			CV%
				< 10%
Sample type	Serum, Plasma, Other biological fluids, Tissue Homogenate			
Assay type	Sandwich (quantitative)			
Sensitivity	< 0.188 ng/ml			
Range	0.313 ng/ml - 20 ng/ml			
Assay duration	Multiple steps standard assay			
Species reactivity	Reacts with: Human			
Product overview	<p>Biosynthesis of L-glutamate from L-aspartate or L-cysteine. Important regulator of levels of glutamate, the major excitatory neurotransmitter of the vertebrate central nervous system. Acts as a scavenger of glutamate in brain neuroprotection. The aspartate aminotransferase activity is involved in hepatic glucose synthesis during development and in adipocyte glyceroneogenesis. Using L-cysteine as substrate, regulates levels of mercaptopyruvate, an important source of hydrogen sulfide. Mercaptopyruvate is converted into H2S via the action of 3-mercaptopyruvate sulfurtransferase (3MST). Hydrogen sulfide is an important synaptic modulator and neuroprotectant in the brain.</p> <p>This Human Aspartate Aminotransferase ELISA Kit (ab285299, E4319) is a sandwich ELISA assay for the quantitative measurement of human AST in serum, plasma and cell culture supernatants. The density of color is proportional to the amount of human AST captured from the samples.</p>			
Notes	This product is manufactured by BioVision, an Abcam company and was previously called E4319			

Aspartate Aminotransferase (AST) (Human) ELISA Kit. E4319-100 is the same size as the 96 test size of ab285299.

Platform Microplate (12 x 8 well strips)

Properties

Storage instructions Store at +4°C. Please refer to protocols.

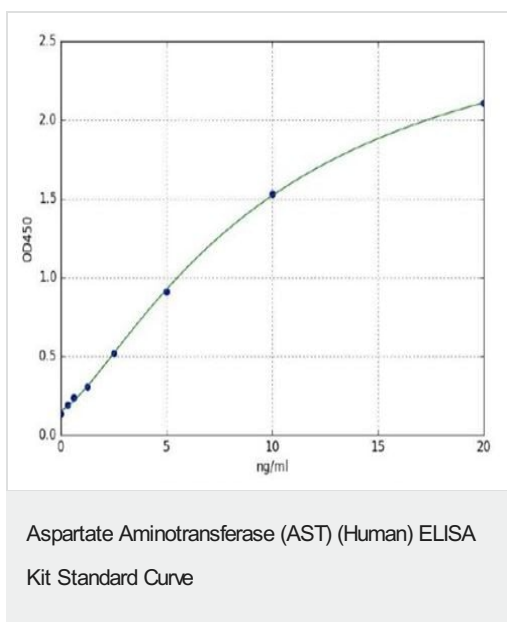
Components	96 tests
25X Wash buffer	1 x 30ml
Antibody dilution buffer	1 x 10ml
Biotin/Detection antibody (Concentrated)	1 x 120µl
HRP-Streptavidin Conjugate (SABC)	1 x 120µl
Lyophilized Standard	2 vials
Micro ELISA Plate	1 unit
Plate sealers	5 units
SABC dilution buffer	1 x 10ml
Sample and Standard Dilution Buffer	1 x 20ml
Stop Solution	1 x 10ml
TMB Substrate	1 x 10ml

Function Plays a key role in amino acid metabolism.

Sequence similarities Belongs to the class-I pyridoxal-phosphate-dependent aminotransferase family.

Cellular localization Cytoplasm.

Images



Typical Standard Curve: These standard curves are for demonstration only. A standard curve must be run with each assay.

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