

Product datasheet

Human CD44var ELISA Kit (v6) ab119519

[1 Image](#)

Overview

Product name Human CD44var ELISA Kit (v6)

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall	8			3%

Inter-assay

Sample	n	Mean	SD	CV%
Overall	8			4.2%

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity 0.05 ng/ml

Range 0.32 ng/ml - 10 ng/ml

Recovery 106 %

Sample specific recovery

Sample type	Average %	Range
Serum	106	89% - 115%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview

Abcam's CD44var (v6) Human in vitro ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for accurate quantitative measurement of Human CD44var (v6) concentrations in Cell culture supernatant, serum and plasma (EDTA, citrate, heparin), and urine.

CD44var (v6) specific antibodies have been precoated onto a 96-well plate. Standards and test samples are added to the wells and along with a CD44var (v6) HRP-conjugated detection antibody and the microplate is then incubated at room temperature. After the removal of unbound proteins by washing, TMB is added and catalyzed by HRP to produce a blue color product that

changes to yellow after addition of an acidic stop solution. The density of yellow coloration is directly proportional to the CD44var (v6) amount of sample captured in plate.

Tested applications

Suitable for: Sandwich ELISA

Platform

Microplate

Properties

Storage instructions

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

Components	1 x 96 tests
20X Assay Buffer Concentrate	1 x 5ml
20X Wash Buffer Concentrate	1 x 50ml
Adhesive Films	2 units
HRP-Conjugate anti-Human CD44var (v6) monoclonal antibody	2 x 10µl
Human CD44var (v6) Standard Lyophilized 20 ng/mL upon reconstitution)	2 vials
Microplate coated with anti CD44var (v6) monoclonal antibody (12 x 8 wells)	1 unit
Sample Diluent	1 x 50ml
Stop Solution (1M Phosphoric acid)	1 x 15ml
TMB Substrate Solution	1 x 15ml

Function

Receptor for hyaluronic acid (HA). Mediates cell-cell and cell-matrix interactions through its affinity for HA, and possibly also through its affinity for other ligands such as osteopontin, collagens, and matrix metalloproteinases (MMPs). Adhesion with HA plays an important role in cell migration, tumor growth and progression. Also involved in lymphocyte activation, recirculation and homing, and in hematopoiesis. Altered expression or dysfunction causes numerous pathogenic phenotypes. Great protein heterogeneity due to numerous alternative splicing and post-translational modification events.

Tissue specificity

Isoform 10 (epithelial isoform) is expressed by cells of epithelium and highly expressed by carcinomas. Expression is repressed in neuroblastoma cells.

Sequence similarities

Contains 1 Link domain.

Domain

The lectin-like LINK domain is responsible for hyaluronan binding.

Post-translational modifications

Proteolytically cleaved in the extracellular matrix by specific proteinases (possibly MMPs) in several cell lines and tumors.

N-glycosylated.

O-glycosylated; contains more-or-less-sulfated chondroitin sulfate glycans, whose number may affect the accessibility of specific proteinases to their cleavage site(s).

Phosphorylated; activation of PKC results in the dephosphorylation of Ser-706 (constitutive phosphorylation site), and the phosphorylation of Ser-672.

Cellular localization

Membrane.

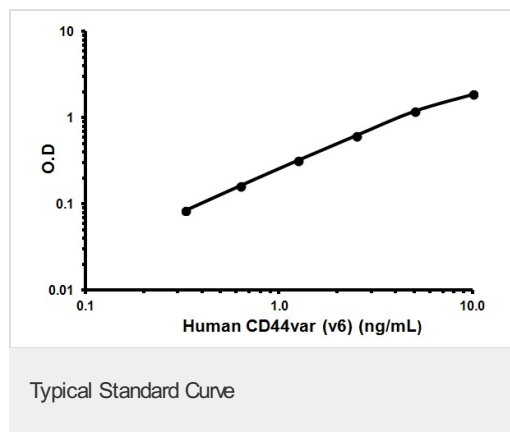
Applications

Our [Abpromise guarantee](#) covers the use of **ab119519** 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
Sandwich ELISA		Use at an assay dependent concentration.

Images



Representative Standard Curve using ab119519.

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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 <http://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