

## 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.

---

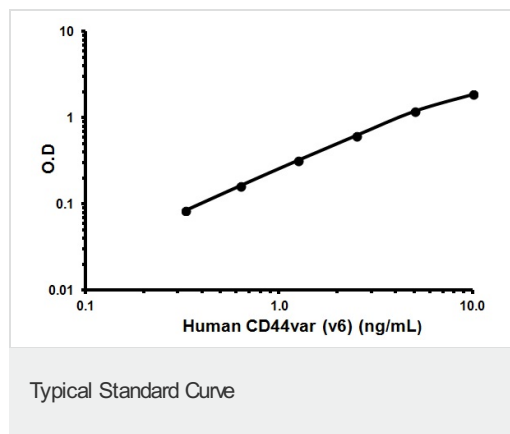
## Applications

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