

### Bradykinin ELISA kit ab136936

★★★★☆ [2 Abreviews](#) [6 References](#) [1 Image](#)

#### Overview

**Product name** Bradykinin ELISA kit

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
Buffer	20	695.4pg/ml		4.6%
Buffer	20	208.7pg/ml		6.2%
Buffer	20	73.7pg/ml		9.9%

Inter-assay

Sample	n	Mean	SD	CV%
Buffer		700.4pg/ml		= 15%
Buffer		209.3pg/ml		= 10.3%
Buffer		66.1pg/ml		= 11.9%

**Sample type** Urine, Serum, Plasma

**Assay type** Competitive

**Sensitivity** 24.8 pg/ml

**Range** 11.7 pg/ml - 30000 pg/ml

**Recovery**

Sample specific recovery

Sample type	Average %	Range
Urine	109	10pg/ml - 2000pg/ml
Serum	113	100pg/ml - 2000pg/ml
Plasma	110	100pg/ml - 20000pg/ml

<b>Assay time</b>	3h 00m
<b>Assay duration</b>	Multiple steps standard assay
<b>Species reactivity</b>	<b>Reacts with:</b> Species independent
<b>Product overview</b>	Bradykinin ELISA kit is a competitive Enzyme-Linked Immunosorbent Assay designed for the accurate quantitative measurement of Bradykinin in plasma, serum and urine.

A goat anti-Rabbit IgG antibody has been precoated onto 96-well plates. Standards or test samples are added to the wells, along with a solution of Bradykinin conjugated to biotin, followed by a solution of polyclonal antibody to Bradykinin. The plate is washed to remove unbound reagents. A solution of streptavidin-HRP conjugate is then added. After further incubation the excess reagents are washed away and TMB substrate is added, which is catalyzed by HRP to generate a yellow color. A stop solution changes this color from yellow to blue, and the intensity of this blue coloration is inversely proportional to the amount of Bradykinin captured in the plate.

**Platform** Microplate

## Properties

**Storage instructions** Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer Concentrate	1 x 27ml
Assay Buffer 16	1 x 55ml
Bradykinin Antibody	1 x 5ml
Bradykinin Conjugate	1 x 5ml
Bradykinin Standard	2 vials
Goat anti-rabbit IgG Microplate (12 x 8 wells)	1 unit
HRP- Streptavidin Conjugate	1 x 12.5µg
Plate Sealer	2 units
Stop Solution 2	1 x 10ml
TMB Substrate	2 x 10ml

## Function

(1) Kininogens are inhibitors of thiol proteases; (2) HMW-kininogen plays an important role in blood coagulation by helping to position optimally prekallikrein and factor XI next to factor XII; (3) HMW-kininogen inhibits the thrombin- and plasmin-induced aggregation of thrombocytes; (4) the active peptide bradykinin that is released from HMW-kininogen shows a variety of physiological effects: (4A) influence in smooth muscle contraction, (4B) induction of hypotension, (4C) natriuresis and diuresis, (4D) decrease in blood glucose level, (4E) it is a mediator of inflammation and causes (4E1) increase in vascular permeability, (4E2) stimulation of nociceptors (4E3) release of other mediators of inflammation (e.g. prostaglandins), (4F) it has a cardioprotective effect (directly via bradykinin action, indirectly via endothelium-derived relaxing

factor action); (5) LMW-kininogen inhibits the aggregation of thrombocytes; (6) LMW-kininogen is in contrast to HMW-kininogen not involved in blood clotting.

**Tissue specificity**

Secreted in plasma. T-kinin is detected in malignant ovarian, colon and breast carcinomas, but not in benign tumors.

**Involvement in disease**

Defects in KNG1 are the cause of high molecular weight kininogen deficiency (HMWK deficiency) [MIM:228960]. HMWK deficiency is an autosomal recessive coagulation defect. Patients with HWMK deficiency do not have a hemorrhagic tendency, but they exhibit abnormal surface-mediated activation of fibrinolysis.

**Sequence similarities**

Contains 3 cystatin domains.

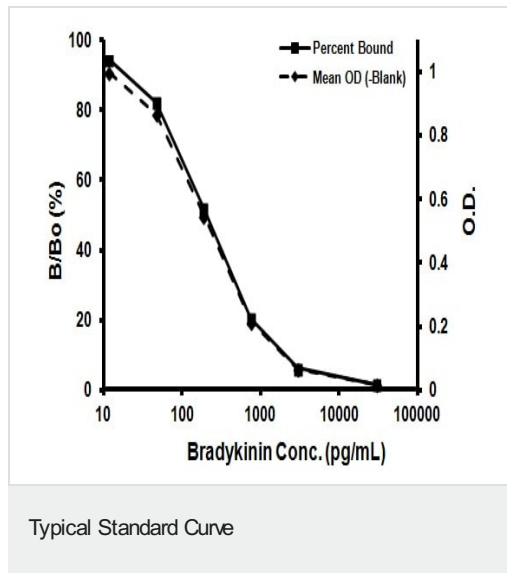
**Post-translational modifications**

Bradykinin is released from kininogen by plasma kallikrein.  
Hydroxylation of Pro-383 occurs prior to the release of bradykinin.  
Phosphorylation sites are present in the extracellular medium.  
N- and O-glycosylated. O-glycosylated with core 1 or possibly core 8 glycans.

**Cellular localization**

Secreted > extracellular space.

**Images**



Representative standard curve using ab136936.

Typical Standard Curve

**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