

Bradykinin ELISA kit

Sensitive and reliable ELISA for measuring bradykinin in plasma, serum and urine samples.

The Bradykinin EIA kit is a colorimetric, competitive immunoassay kit with results in 3 hours.

Bradykinin was discovered in 1949 as a substance generated from a globulin precursor in plasma by the action of proteases. Its name indicates that it causes a slow movement of the gut. As early as 1909 it was noted that substances found in urine, which were later identified as kinins, have hypotensive actions. Kinins are effectors of vasodilation, vascular permeability, NO release and arachidonic acid mobilization. They are important regulators of blood pressure, kidney function and heart function, and they are also involved in inflammation. Bradykinin is generated from the blood globulin kininogen HK, by the action of the kallikrein system in blood (related to the blood clotting cascade) but can also be generated in other tissues and organs. Besides kallikrein, other proteases such as plasmin may also release bradykinin. Several peptidases can degrade kinins, including Angiotensin Converting Enzyme (ACE), a metalloproteinase which converts Angiotensin I to Angiotensin II and destroys bradykinin. Plasma bradykinin is rapidly degraded to a smaller stable peptide (BK1-5) form.

Citations: 22

[View Online »](#)

Ordering Information

[Order Online »](#)

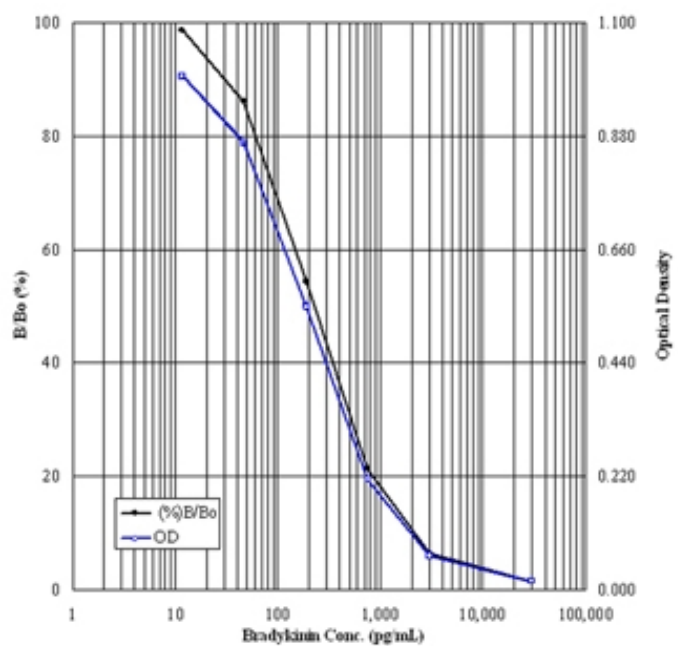
ADI-900-206

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Sensitive, reliably measure as little as 24.8 pg/ml of Bradykinin
- Rapid results in just 3 hours from up to 41 samples in duplicate
- Broad dynamic range to easily quantify high and low levels of Bradykinin
- Easy-to-use liquid color-coded reagents reduce errors



Handling & Storage

| | |
|--------------------|---|
| Use/Stability | Store all components at +4°C, except Standard at -20°C. |
| Short Term Storage | -20°C |
| Long Term Storage | -20°C |
| Shipping | Blue Ice |

Regulatory Status

RUO - Research Use Only

Product Details

| | |
|--------------------|---|
| Application | Colorimetric detection, ELISA |
| Application Notes | For the quantitative determination of species independent Bradykinin in plasma, serum and urine samples. |
| Assay Time | 3 hours |
| Compatibility | This product is compatible with the Absorbance 96 Plate Reader . |
| Contents | Microtiter Plate, Conjugate, Antibody, Assay Buffer 16, Wash Buffer Concentrate, Standard, TMB Substrate, Stop Solution 2, Streptavidin-HRP |
| Crossreactivity | Bradykinin (100%) Lys-Bradykinin (Kallidin) (100%) and Les-des-Arg9-Bradykinin (<1%), BK1-5 stable degradion product (<0.1%) |
| Quantity | 1 x 96-well plate |
| Sensitivity | 24.8 pg/ml (range 11.7-30,000 pg/ml) |
| Species Reactivity | Species independent |
| UniProt ID | P01042 (human) |
| Wavelength | 450 nm |



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)