

BNP Fragment ELISA kit

Sensitive BNP fragment ELISA kit enabling detection of BNP fragment in serum and plasma.

The BNP Fragment ELISA kit is a colorimetric competitive enzyme immunoassay kit.

Brain natriuretic peptide (BNP) is secreted by the ventricles of the heart in response to excessive stretching of cardiomyocytes. The physiologic actions of BNP include decrease in systemic vascular resistance and central venous pressure and an increase in natriuresis. BNP measurements are used to aid in the diagnosis and assessment of severity of heart failure.

Ordering Information

[Order Online »](#)

ENZ-KIT152-0001	96 wells
-----------------	----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	All reagents of the kit are stable at 4°C (2-8°C) until the expiry date.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Brain natriuretic peptide
Application	Colorimetric detection, ELISA
Application Notes	For the quantitative determination of human BNP fragment in serum and plasma.
Assay Time	O/N plus 20 mins
Compatibility	This product is compatible with the Absorbance 96 Plate Reader .
Sensitivity	171 pmol/l (0-6400 pmol/l)
Species Reactivity	Human
UniProt ID	P16860
Wavelength	450 nm

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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