

NT-proCNP ELISA kit

Sensitive NT-proCNP ELISA kit enabling detection of NT-proCNP in serum and plasma in just 4 hours.

The NT-proCNP ELISA kit is a colorimetric immunometric enzyme immunoassay kit with results in 4 hours.

C-type natriuretic peptide (CNP) is a hormone which plays a role in endochondral ossification through regulation of cartilaginous growth plate chondrocytes proliferation and differentiation. CNP may also be vasoactive and natriuretic.

Ordering Information

[Order Online »](#)

ENZ-KIT154A-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 C-type natriuretic peptide

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of human NT-proCNP in serum, EDTA plasma, heparin plasma, and citrate plasma.

Assay Time 4 hours

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

Sensitivity 0.7 pmol/l (0-128 pmol/l)

Species Reactivity Human

UniProt ID P23582

Wavelength 450 nm

Last modified: May 29, 2024



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)