

Procollagen type I (CT) (human) monoclonal antibody (4)

Type I collagen is present in connective tissue and bone where it makes up about 90% of the organic matrix. Type I collagen is produced from type-I procollagen by the proteolytic cleavage of N- and C-telopeptides, which are released into the circulation. The propeptides can be measured in serum as markers of the production of type-I collagen. Elevated values of these propeptides are found in serum from patients with various metabolic or malignant bone diseases. Procollagen type I C-terminal propeptide (PICP) has a molecular mass of approximately 100kDa. Its normal concentration in serum depends on age and sex.

This antibody is covered by our [Worry-Free Guarantee](#).

Ordering Information

[Order Online »](#)

BPD-BTE-002-04-1	1mg
BPD-BTE-002-04-04	400µg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Type I collagen, Procollagen type I C-terminal propeptide, PICP

Application ELISA

Clone 4

Formulation Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.

GenBank ID 1277

Host Mouse

Immunogen Purified human PICP from second trimester amniotic fluid.

Isotype IgG1κ

Purity Detail Protein A/G-purified.

Source From culture supernatant.

Species Reactivity Human

Specificity Specific for native PICP from amniotic fluid.

UniProt ID P02452

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#).



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)