

Procollagen type I (CT) (human) monoclonal antibody (8)

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-002-08-1	1mg
BPD-BTE-002-08-04	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, IP
Clone	8
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 for native PICP from amniotic fluid.
UniProt ID	P02452

Worry-free Guarantee

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