

# 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](#).

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](#)



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)

Last modified: May 29, 2024