# 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

Purified human PICP from second trimester amniotic fluid. **Immunogen** 

lgG1κ Isotype

**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



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 usa@enzolifesciences.com eu@enzolifesciences.com Last modified: May 29, 2024

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

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

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

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