

# Procollagen type IIA (NT) (human) monoclonal antibody (2)

Collagen type II is the most abundant protein of cartilage matrix. Two isoforms exist, and collagen type IIA is involved in early development of cartilage in embryos. The collagen is produced from type II procollagen when two teloterminal propeptides are digested from the precursor and released into the circulation. Procollagen type II N-terminal propeptide (PIIANP) is considered a marker of cartilage synthesis, and the normal concentration in serum is approximately 30 ng/mL.

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

## Ordering Information

[Order Online »](#)

BPD-BTE-003-02-04	400µg
-------------------	-------

## Manuals, SDS & CofA

[View Online »](#)

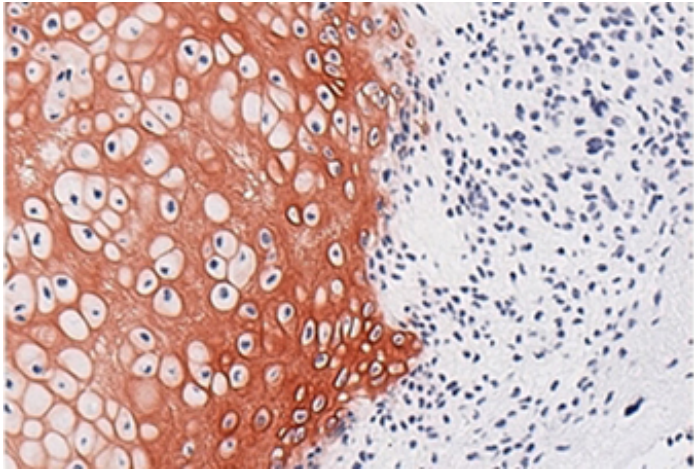


Figure 2: Immunohistochemical analysis of formalin-fixed, paraffin-embedded biopsy from a chondroblastoma, showing specific staining of tumor cells using Procollagen type IIA (NT) (human), mAb and developed with peroxidase-labeled goat anti-mouse Ig (blue color).

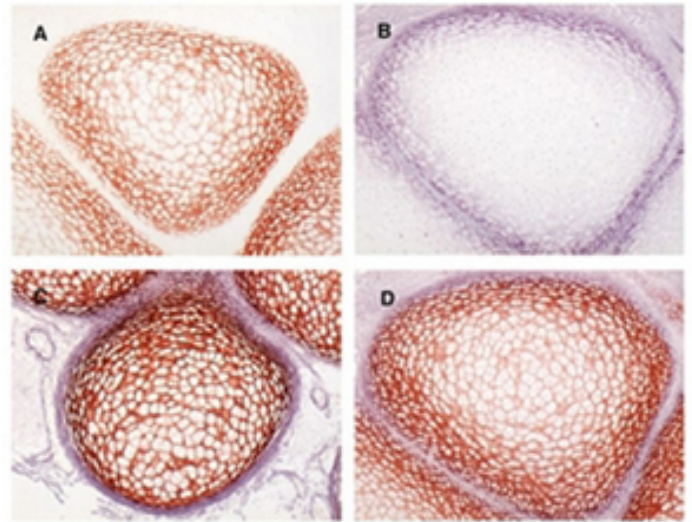


Figure 1: Immunohistochemical analysis of formalin-fixed, paraffin-embedded fetal bone anlage in the hand (around week 10) stained using Procollagen type IIA (NT).

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	Collagen type II, Procollagen type II N-terminal propeptide, PIIANP
Application	ELISA, IHC (PS)
Clone	2
Formulation	Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.
GenBank ID	1280
Host	Mouse
Immunogen	Recombinant PIIANP.
Isotype	IgG1
Purity Detail	Protein A/G-purified.
Source	From culture supernatant.
Species Reactivity	Human
Specificity	Epitope is located at the amino-terminal part of procollagen type IIA.
UniProt ID	P02458
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>

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)