

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-003-02-1	1mg
BPD-BTE-003-02-04	400µg

Manuals, SDS & CofA

[View Online »](#)

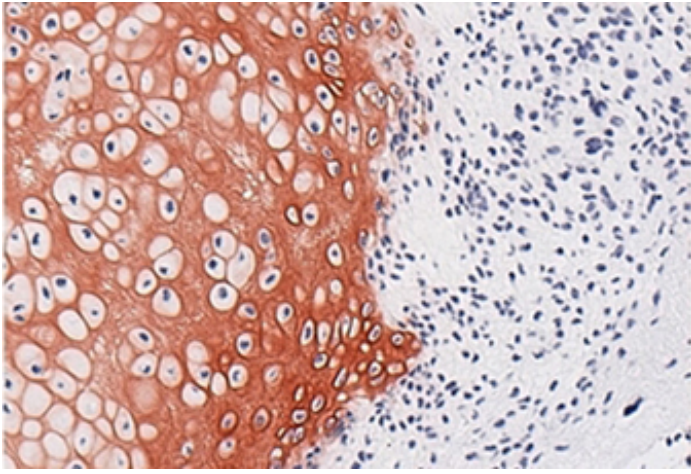


Figure 2: Immunohistochemical analysis of formalin-fixed, paraffin-embedded biopsy from a chondroblastoma, showing specific staining of tumor cells using Procollagen type IIA (NT) (human), mAb and developed with peroxidase-labeled goat anti-mouse Ig (blue color).

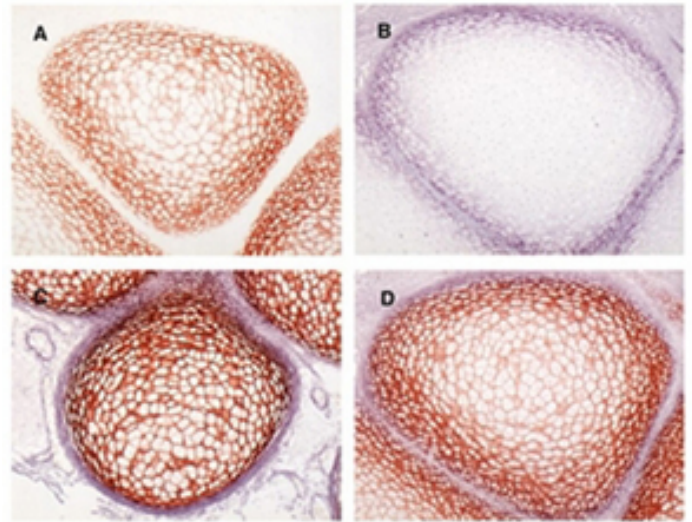


Figure 1: Immunohistochemical analysis of formalin-fixed, paraffin-embedded fetal bone anlage in the hand (around week 10) stained using Procollagen type IIA (NT).

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 [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)