

Collagen Type V monoclonal antibody (1E2-E4/Col5)

Type V collagen is a minor component of the connective tissue, although it is present in many different types of connective tissue. Patients with defects in the type V collagen (Ehlers-Danlos syndrome) have weakened connective tissue characterized by hyperstretchable joints and fragile, easily bruisable skin.

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

Ordering Information

[Order Online »](#)

BPD-CSI-006-01-1	1mg
BPD-CSI-006-01-04	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 COL5

Application ELISA, IHC (FS), IHC (PS), WB

Clone 1E2-E4/Col5

Crossreactivity Does not cross-react with mouse, rat, guinea pig or chicken. It has been shown to have no cross-reactivity with type I, III and VI collagens by ELISA and immunoblotting. There is no evidence for cross-reactivity with other connective tissue proteins (laminin, fibronectin, elastin).

Formulation Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.

GenBank ID 1289

Host Mouse

Immunogen Acid-digested pepsin soluble dog type V collagen.

Isotype IgG2ak

Purity Detail Protein A/G-purified.

Source From culture supernatant.

Species Reactivity Bovine, Dog, Human, Kangaroo, Porcine, Rabbit, Sheep

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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-fr@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)