

# 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](#).

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 <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)