

# Collagen type XII monoclonal antibody (378D5)

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

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Application	ELISA, IHC (FS), IP, WB
Application Notes	Not recommended for IHC (PS).
Clone	378D5
Formulation	Liquid. Ascites. Contains no preservatives.
Host	Mouse
Immunogen	Collagen type XII purified from bovine kidney.
Isotype	IgG1
Recommendation Dilutions/Conditions	ELISA (1:500-1:1,000)Immunohistochemistry (frozen sections, 1:500-1:1,000; a mild trypsin- or Protease 27 (Sigma)-digestion (0.05% for 10-15 min at 37°C) or alternatively microwave oven treatment is recommended to unmask the collagens)Immunoprecipitation (1:50-1:100)Western Blot (1:100-1:300)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Bovine, Chicken, Human, Mouse, Rat
UniProt ID	P25508
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)