

Collagen type XII monoclonal antibody (378D5)

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

Ordering Information

[Order Online »](#)

ALX-803-309-R100	100µl
------------------	-------

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 [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)