

Anti-Collagenase (Clostridium histolyticum), polyclonal antibody from sheep

Ordering Information

[Order Online »](#)

| | |
|-------------|-------|
| NB-58050-01 | 100µg |
|-------------|-------|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ELISA, IHC, WB

Formulation In PBS buffer with 0.01 % sodium azide.

Host Sheep

Immunogen Clostridial Collagenase NB.

Isotype IgG

Species Reactivity Clostridium histolyticum

Specificity Recognizes several collagenase isoforms.

Technical Info / Product NB-58050-01 corresponds to supplier's Cat. No. S5805001

Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com