

Caspase-3 polyclonal antibody

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

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

| | |
|------------------|-------|
| ALX-210-806-C100 | 100µg |
|------------------|-------|

Manuals, SDS & CofA

[View Online »](#)



Figure: Western blot analysis of human caspase-3 using PAb to caspase-3 (Prod. No. ALX-210-806) with (lane 1) and without (lane 2) blocking peptide.

Handling & Storage

| | |
|-------------------|---|
| Handling | Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C. |
| Long Term Storage | -20°C |
| Shipping | Blue Ice |

Regulatory Status

RUO - Research Use Only

Product Details

| | |
|-------------------------------------|--|
| Application | WB |
| Application Notes | Detects bands of ~32kDa (procaspase-3) and ~20kDa (cleavage product) by Western blot. |
| Formulation | Liquid. In PBS, pH 7.2, containing 0.5% BSA, 5mM EDTA, 0.03% Proclin and 30% glycerol. |
| Host | Rabbit |
| Immunogen | Synthetic peptide corresponding to aa 163-175 of human caspase-3. |
| Purity Detail | Epitope-affinity purified. |
| Recommendation Dilutions/Conditions | Western Blot (0.5-4µg/ml) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application. |
| Species Reactivity | Human, Mouse, Rat |
| UniProt ID | P42574 |
| Worry-free Guarantee | This antibody is covered by our Worry-Free Guarantee |

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)