

Procaspase-9 (mouse), (recombinant)

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-201-165-C050	50µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability Stability at -20°C: 1-2 weeks.

Handling Avoid freeze/thaw cycles.

Short Term Storage -20°C

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Concentration 0.57mg/ml

Formulation Liquid. In 20mM TRIS-HCl, pH 8.5.

Purity ≥90% (SDS-PAGE)

Source Produced in *E. coli*. Requires activation.

UniProt ID Q8C3Q9

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-
fr@enzolifesciences.com](mailto:info-fr@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)