

Caspase-9 (human), (recombinant)

Caspase-9 is the initiator caspase of the mitochondrial/Apaf-1 apoptosis induction pathway.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE173-5000	5000U
----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Activity 100 U/μl

Alternative Name ICE-LAP-6, Mch6

Formulation Liquid. In 50mM HEPES, pH 7.4, 100mM NaCl, 0.5% CHAPS, 1mM EDTA, 10% glycerol, 10mM DTT.

MW 20 + 12 kDa

Purity ≥90% (SDS-PAGE)

Source Produced in *E. coli*. C-terminal Human caspase-9 (aa 131-416).

Specific Activity One U=1 pmol/min, at 30°C using Ac-LEHD-pNA (2mM, Prod. No. ALX-260-081) as substrate when measuring 100 U diluted in assay buffer consisting of 100mM MES pH 6.5, 10% PEG 8000, 0.1% CHAPS, 10 MM DTT.

UniProt ID P55211

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)