

Caspase-4 (human), (recombinant)

Caspase-4 is localized primarily to the endoplasmic reticulum (ER) and is required for mediation of ER-stress induced apoptosis.

Citations: 9

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE176-5000	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	ICH-2, Protease TX, ICereI-II
Formulation	Liquid. In 50mM HEPES, pH 7.4, 100mM NaCl, 0.5% CHAPS, 1mM EDTA, 10% glycerol, 10mM DTT.
MW	19 + 10 kDa
Purity	≥90% (SDS-PAGE)
Source	Produced in <i>E. coli</i> .
Specific Activity	One U=1 pmol/min at 30°C using Ac-LEHD-pNA (200 mM; Prod. No. ALX-260-081) as substrate.
UniProt ID	P49662

Last modified: May 29, 2024



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