

Caspase-5 (human), (recombinant)

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE171-5000	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-3, Protease TY, ICErel-III
Formulation	Liquid. In 50mM HEPES, pH 7.4, 100mM NaCl, 0.5% CHAPS, 1mM EDTA, 10% glycerol, 10mM DTT.
MW	22 + 10 kDa
Purity	≥80% (SDS-PAGE)
Source	Produced in <i>E. coli</i>
Specific Activity	One U=1 pmol/min at 30°C using Ac-LEHD-pNA (200 μM; Prod. No. ALX-260-081) as substrate.
UniProt ID	P51878

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)