

Caspase-11 (mouse), (recombinant)

A member of the caspase-1 subfamily, caspase-11 is the possible homolog of human caspase-4. Its expression is highly inducible, in various mouse tissues, by LPS injection.

Citations: 5

[View Online »](#)

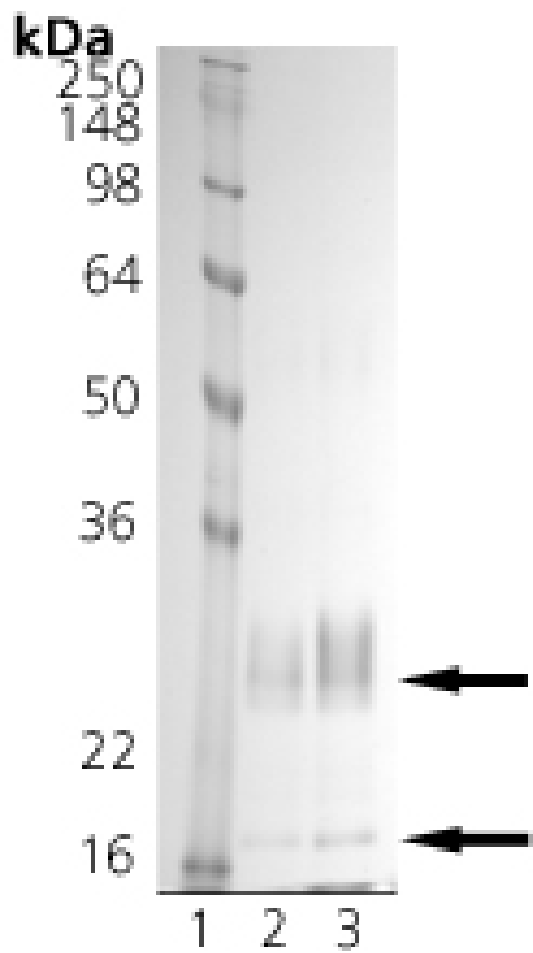
Ordering Information

[Order Online »](#)

BML-SE155-5000	5000U
----------------	-------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis: Lane 1: MW Marker, Lane 2: 1 μ g, Lane 2: 2 μ g Caspase-11.

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Activity	100 U/μl
Formulation	Liquid. In 50mM HEPES, pH 7.4, 100mM NaCl, 0.5% CHAPS, 1 mM EDTA, 10% glycerol, 10mM DTT.
MW	23 + 10 kDa subunits
Purity	≥95% (SDS-PAGE)
Source	Produced in <i>E. coli</i> . Mouse caspase-11 (aa 81-373).
Specific Activity	One unit hydrolyzes 1 pmol/min of Ac-WEHD-pNA (500 μM, Prod. No. ALX-260-082) at 30°C.
UniProt ID	P70343

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)