

Caspase-10 (human), (recombinant)

Caspase-10 is an initiator caspase, highly homologous to caspase-8, a relationship that extends to the death effector domains (DEDs) of their proforms. Caspase-10 was reported to activate early in TRAIL-induced apoptosis and was found in association with the DR4 (TRAIL-R1) receptor complex in a thyroid carcinoma line.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE174-5000	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 FLICE2, Mch4

Formulation Liquid. In 50mM HEPES, pH 7.4, 100mM NaCl, 0.1% CHAPS, 1mM EDTA, 10% glycerol, 10mM DTT.

MW 17 + 12 kDa

Purity ≥95% (SDS-PAGE)

Source Produced in *E. coli*.

Specific Activity One U=1pmol/min using Ac-IETD-pNA (200μM; Prod. No. ALX-260-045) as substrate at 30°C.

UniProt ID Q92851



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)