Caspase-8 (human), (recombinant)

Highly active caspase-8 for the study of enzyme regulation and kinetics

Caspase-8 is an initiator caspase activated in the death-inducing signaling complexes (DISCs) of the TNF receptor superfamily (e.g. TRAIL, FasL, TNF α receptors).

Citations: 8

View Online »

Ordering Information

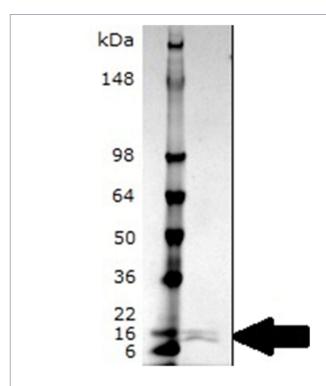
Order Online »

BML-SE172-5000

5000U

Manuals, SDS & CofA

View Online »



SDS PAGE Analysis: Lane 1: MW Marker, Lane 2: 1 μ g (Prod. No. BML-SE172).

Handling & Storage

Use/Stability After initial defrost, aliquot product into individual tubes and refreeze the remaining,

unused enzyme quickly by snap-freezing in a dry/ice ethanol bath or liquid nitrogen, if possible. Avoid repeated freeze/defrost cycles.NOTE: When stored under the above conditions, this enzyme is stable at the concentration supplied, in its current storage buffer. Procedures such as dilution of the enzyme followed by refreezing, could lead to

loss of activity.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Activity 100 U/µI

Alternative Name FLICE, Mch5, MACH

Application NotesUseful tool to study enzyme regulation and kinetics, cleave

target substrates, screen for inhibitors.

Formulation Liquid. In 50mM HEPES, pH 7.4, 100mM sodium chloride,

0.5% CHAPS, 1mM EDTA, 10% glycerol and 10mM DTT.

MW 18 + 11 kDa

Purity ≥90% (SDS-PAGE)

Source Produced in E. coli

Specific Activity One U=1 pmol/min using Ac-IETD-pNA (200mM; Prod.

No. ALX-260-045) as substrate, at 30°C.

UniProt ID Q14790

Last modified: May 29, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com