Caspase-6 (human), (recombinant)

Highly, active caspase-6 for the study of enzyme kinetics or screening of inhibitors

Effector caspase responsible for the cleavage of nuclear lamins in apoptosis. Caspase-6 cleavages of various other protein targets have been implicated in the pathogenesis of neurodegenerative diseases, including Alzheimer's disease (b-amyloid precursor protein), Huntington's disease (huntingtin).

Citations: 10

View Online »

Ordering Information

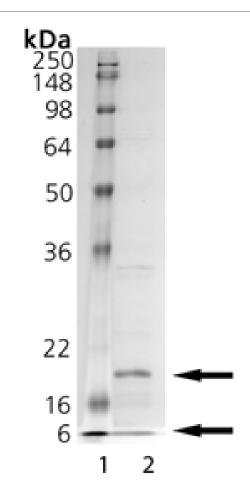
Order Online »

BML-SE170-5000

5000U

Manuals, SDS & CofA

View Online »



SDS-PAGE Analysis – Lane 1: MW Marker, Lane 2: 1 µg Caspase-6.

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.

Handling Avoid freeze/thaw cycles.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Activity 100 U/μl

Alternative Name Mch2

Application Notes Useful tool to study enzyme regulation and kinetics, cleave

target substrates, screen for inhibitors.

Formulation Liquid. In 50mM HEPES, pH 7.4, containing 100mM

sodium chloride, 0.5% CHAPS, 1mM EDTA, 10% glycerol

and 10mM DTT.

MW 18 + 11 kDa subunits

Purity ≥80% (SDS-PAGE)

Source Produced in *E. coli*.

Specific Activity One U=1 pmol/min at 30°C using Ac-VEID-pNA (200mM;

Prod. No. BML-P437) as substrate.

UniProt ID P55212

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