Caspase-8 (human), (recombinant) (active)

Caspase-8 is a member of the interleukin-1 β converting enzyme (ICE) family of cysteine proteases. Same as other caspases, caspase-7 also exists in cells as an inactive proenzyme. During apoptosis procaspase-8 is processed at aspartate residues by self-proteolysis and/or cleavage by another caspase. The processed form of caspase-8 consists of large and small subunits which associate to form the active enzyme.

Citations: 5

View Online »

Ordering Information

Order Online »

ALX-201-062-U100

100U

Manuals, SDS & CofA

View Online »

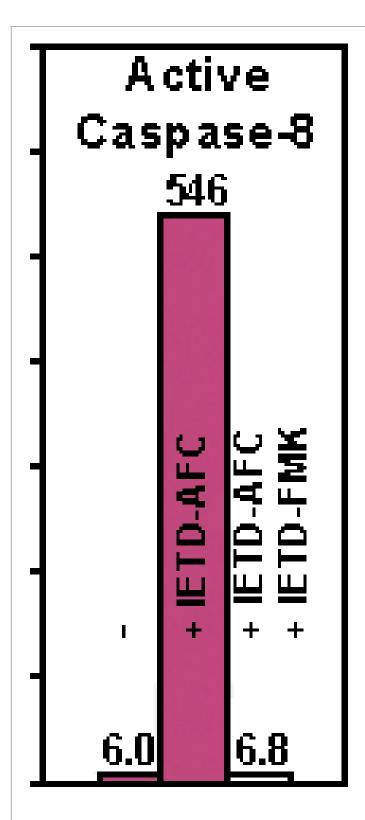


Figure: Active human caspase was expressed in *E. coli* and purified. The activity of recombinant caspase-8 was determined by cleaving AFC conjugates of IETD. The cleavage activity was effectively inhibited by the corresponding peptide inhibitor as indicated.

Handling & Storage

Handling Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name FLICE, Mch5, MACH

Application NotesUseful in screening caspase inhibitors, studying enzyme

regulation, determining specificity of caspase substrates or

as positive control in caspase activity assays. We recommend using 1U per assay for analyzing caspase

activity.

Formulation Lyophilized.

Purity ≥95% (SDS-PAGE)

Reconstitution Reconstitute to 1U/µl with PBS containing 15% glycerol.

Source Produced in *E. coli*. Contains an N-terminal His-tag.

Specific Activity 5'000U/mg protein. One unit is defined as the amount of

enzyme that cleaves 1nmol of the caspase substrate IETD-pNA per hour at 37°C in a reaction solution

containing 50mM HEPES, pH 7.2, 50mM sodium chloride,

0.1% CHAPS, 10mM EDTA, 5% glycerol and 10mM DTT.

UniProt ID Q14790

Last modified: May 29, 2024

