Caspase-4 (human), (recombinant) (active)

Caspase-4 is a member of the caspase family of cysteine proteases. It exists in cells as an inactive pro-enzyme. The pro-enzyme is matured by proteolysis to yield large and small subunits. The active enzyme is a heterotetramer consisting of two large and two small subunits. The expressed caspase-4 spontaneously undergoes auto-processing to yield subunits characteristic of the native enzyme. The active caspase-4 is routinely tested for its ability to enzymatically cleave the substrate WEHD-AFC or WEHD-pNA.

Citations: 2

View Online »

Ordering Information

Order Online »

ALX-201-093A-U025	25U
ALX-201-093A-U100	100U

Manuals, SDS & CofA

View Online »

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 ICH-2, Protease TX, ICErel-II

Application NotesUseful in screening caspase inhibitors, studying enzyme

regulation and kinetics, determining target substrate or as

a positive control in caspase activity assays. We

recommend using 1 unit per assay for analyzing caspase

activity.

Formulation Lyophilized.

Purity ≥95% (SDS-PAGE)

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

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 WEHD-pNA per hour at 37°C in reaction solution containing 50mM HEPES, pH 7.2, 50mM NaCl, 0.1% CHAPS, 10mM EDTA, 5% glycerol and 10mM DTT.

UniProt ID P49662

Last modified: May 29, 2024

