

Instruction Manual

Catalog Number	PK-RP577-1181-100
Description	Caspase-1 (also known as ICE) is a prototypical member of the caspase-family of cysteine proteases. Caspase-1 exists in cells as an inactive 45 kDa proenzyme. The pro-enzyme is matured by proteolysis to yield large (20 kD) and small (10 kD) subunits. The active caspase-1 is a heterotetramer consisting of two large and two small subunits. To date the regulatory mechanism of caspase-1 activation and the role of caspase-1 in apoptosis are poorly understood. In THP-1 cells, a large proportion of the caspase-1 is present in the inactive proenzyme form. The recombinant active mouse caspase-1 was expressed in E. coli. The active caspase-1 preferentially cleaves caspase-1 substrates (e.g., YVAD-AFC or YVAD-pNA) and is routinely tested for its ability to enzymatically cleave these two substrates Ac-YVAD-pNA or Ac-YVAD-AFC.
Quantity	100 units
Specific Activity	>3,500 units/mg
Unit Definition	One unit of the recombinant caspase-1 is the enzyme activity that cleaves 1 nmol of the caspase substrate Ac-YVAD-pNA (pNA: p-nitroaniline) per hour at 37°C in a reaction solution containing 50 mM HEPES, pH 7.2, 50 mM NaCl, 0.1% CHAPS, 10 mM EDTA, 5% Glycerol, and 10 mM DTT.
Formulation	Lyophilized powder
Reconstitution	Reconstitute to 1 unit per µl in PBS containing 15% glycerol.
Storage	The lyophilized caspase-1 is stable for 1 year at -70°C. Following reconstitution in PBS, the enzyme should be aliquoted and immediately stored at -70°C. Avoid multiple freeze/thaw cycles as activity might decrease.
Applications	Active caspase-1 is useful in studying enzyme regulation, determining target substrates, screening caspase inhibitors, or as a positive control in caspase activity assays. We recommend using 1 unit/assay for analyzing caspase activity. For a complete caspase-1 assay protocol, please refer to PromoKine's Caspase-1 Fluorometric or Colorimetric Assay Kits.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.