

## Instruction Manual

<b>Catalog Number</b>	PK-RP577-1089P-50
<b>Description</b>	Caspase-9 is a member of the caspase family of cysteine proteases. Caspase-9 exists in cells as an inactive proenzyme. Proteolytic processing of this inactive proenzyme generates the large and small subunits which, when assembled as a tetramer (a pair of heterodimers), form the active caspase. Cascades of caspase activation have been shown to be the key signal-transducing events in apoptosis. Caspase-9 is considered to be an upstream enzyme that initiates caspase cascade leading to apoptosis. The recombinant Procaspase-9 is produced by expression of a human cDNA in E. coli and further purified by multiple chromatographic steps.
<b>Quantity</b>	50 µg
<b>Unit Definition</b>	NA
<b>Purity</b>	>90% by SDS-PAGE
<b>Formulation</b>	Freeze-dried powder from 50 mM Tris-HCl buffer, pH 8.0.
<b>Storage</b>	Store at -80°C. Aliquot product into individual vials and freeze immediately at -80°C. AVOID FREEZE/THAW CYCLES!
<b>Applications</b>	Investigating mechanisms of the proteolytic processing and activation of procaspase-9.

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