

## Caspase-9

(CASP-9, ICE-like apoptotic protease 6, ICE-LAP6, Apoptotic protease Mch-6, Apoptotic protease-activating factor 3, APAF-3)

- **CAT #:** Caspase9
- **Enzyme:** Caspase-9, Highly active casp-9 from human cDNA, produced in an *E. coli* expression system (Residues Ile 131 to Ser 416)
- **Substrate:** (Z-LEHD)<sub>2</sub>-R110 tetrapeptide
- **References Compound IC50s:**

<u>Compounds</u>	<u>Ac-IETD-CHO</u>	<u>Ac-LEHD-CHO</u>
IC50(M)	9.79E-08	5.72E-08

- **Scientific Information:** [P55211](#)

\* R110 = Rhodamine 110

### Control inhibitor IC50s for Caspase 9

