

## Caspase-7

(CASP-7, ICE-like apoptotic protease 3, ICE-LAP3 Apoptotic protease Mch-3, CMH-1)

- **CAT #:** Caspase7
- **Enzyme:** Caspase-7, from human cDNA produced in an *E. coli* expression system (residues Ala-24 to Gln-303)
- **Substrate:** (Z-DEVD)<sub>2</sub>-R110 tetrapeptide
- **References Compound IC50s:**

<u>Compounds</u>	<u>Ac-VEID-CHO</u>	<u>Ac-IETD-CHO</u>	<u>Ac-DEVD-CHO</u>
<u>IC50(M)</u>	1.14E-07	1.07E-06	3.35E-09

- **Scientific Information:** [P55210](#)

\* R110 = Rhodamine 110

### Control inhibitor IC50s for Caspase 7

