

## Caspase-6

(CASP-6, Apoptotic protease Mch-2)

- **CAT #:** Caspase6
- **Enzyme:** Caspase-6, from human cDNA produced in an *E. coli* expression system (residues Ala-24 to Asn-293)
- **Substrate:** (Z-DEVD)<sub>2</sub>-R110 tetrapeptide
- **References Compound IC50s:**

<u>Compounds</u>	<u>Ac-IETD-CHO</u>	<u>Ac-DEVD-CHO</u>
<b>IC50(M)</b>	3.11E-08	2.82E-08

- **Scientific Information:** [P55212](#)

\* R110 = Rhodamine 110

### Control inhibitor IC50s for Caspase 6

