

Caspase-3

(CASP-3, Apopain, Cysteine protease CPP32, Yama protein, CPP-32 SREBP cleavage activity 1, SCA-1)

- **CAT #:** Caspase3
- **Enzyme:** Caspase-3, from human cDNA produced in an *E. coli* expression system
- **Substrate:** (Z-DEVD)₂-R110 tetrapeptide
- **References Compound IC50s:**

<u>Compounds</u>	<u>Ac-VEID-CHO</u>	<u>Ac-IETD-CHO</u>
IC50(M)	2.08E-08	5.39E-07

Scientific Information: [P42574](#)

*R110 = Rhodamine 110

Control inhibitor IC50s for Caspase 3

