

Caspase-10

(CASP-10, ICE-like apoptotic protease 4, Apoptotic protease Mch-4, FAS-associated death domain protein interleukin-1B-converting enzyme 2, FLICE2)

- **CAT #:** Caspase10
- **Enzyme:** Caspase-10, Highly active casp-10 from human cDNA, produced in an *E. coli* expression system (Residues Val 220 to Ile 479)
- **Substrate:** (Z-LEHD)₂-R110 tetrapeptide
- **References Compound IC50s:**

<u>Compounds</u>	<u>Ac-IETD-CHO</u>	<u>Ac-VEID-CHO</u>
IC50(M)	1.73E-06	5.34E-07

- **Scientific Information:** [Q92851](#)

* R110 = Rhodamine 110

Control inhibitor IC50s for Caspase 10

