

Caspase-8

(CASP-8, ICE-like apoptotic protease 5, MORT1-associated CED-3 homolog, MACH, FADD-homologous ICE/CED-3-like protease, FADD-like ICE, FLICE, Apoptotic cysteine protease, Apoptotic protease Mch-5, CAP4)

- **CAT #:** Caspase8
- **Enzyme:** Caspase-8, Highly active casp-8 from human cDNA, produced in an *E. coli* expression system (Residues Ser 217 to Asp 479)
- **Substrate:** (Z-LEHD)₂-R110 tetrapeptide
- **References Compound IC50s:**

<u>Compounds</u>	<u>Ac-YVAD-CHO</u>	<u>Ac-DEVD-CHO</u>
IC50(M)	3.94E-08	1.46E-09

- **Scientific Information:** [Q14790](#)

* R110 = Rhodamine 110

Control inhibitor IC50s for Caspase 8

