

Caspase-4

(CASP-4, ICH-2 protease, TX protease, ICE(rel)-II)

- **CAT #:** Caspase4
- **Enzyme:** Caspase-4, Highly active casp-4 from human cDNA, produced in an *E. coli* expression system (Residues Ala 105 to Asn 377)
- **Substrate:** (Z-LEHD)₂-R110 tetrapeptide
- **References Compound IC50s:**

<u>Compounds</u>	<u>Ac-VEID-CHO</u>	<u>Ac-YVAD-CHO</u>	<u>Ac-LEHD-CHO</u>
IC50(M)	2.25E-06	2.37E-06	9.71E-07

- **Scientific Information:** [P49662](#)

* R110 = Rhodamine 110

Control inhibitor IC50s for Caspase 4

