

## Caspase-1

(CASP-1, Interleukin-1 beta convertase, IL-1BC, IL-1 beta-converting enzyme, ICE, Interleukin-1 beta-converting enzyme, p45)

- **CAT #:** Caspase1
- **Enzyme:** Caspase-1, Highly active caspase-1 (ICE) from human cDNA, produced in an *E. coli* expression system (residues 120-404, Asp-381 to Glu change)
- **Substrate:** (Z-LEHD)<sub>2</sub>-R110 tetrapeptide
- **References Compound IC50s:**

<u>Compounds</u>	<u>Ac-VEID-CHO</u>	<u>Ac-YVAD-CHO</u>	<u>Ac-DEVD-CHO</u>
IC50(M)	7.90E-08	2.83E-09	1.02E-06

- **Scientific Information:** [P29466](#)

\* R110 = Rhodamine 110

### Control inhibitor IC50s for Caspase 1

