

## Caspase-14

(CARD-containing MAGUK protein 2, Carma 2)

- **CAT #:** Caspase14
- **Enzyme:** Caspase-14, Highly active casp-14, produced by treatment of procaspase-14 with granzyme B
- **Substrate:** (Z-LEHD)<sub>2</sub>-R110 tetrapeptide
- **References Compound IC50s:**

<u>Compounds</u> IC50(M)	<u>Ac-LEHD-CHO</u> 7.90E-07	<u>Ac-VEID-CHO</u> 2.93E-06	<u>Ac-YVAD-CHO</u> 9.89E-08
-----------------------------	--------------------------------	--------------------------------	--------------------------------

- **Scientific Information:** [Q9BXL6](#)

\* R110 = Rhodamine 110

### Control inhibitor IC50s for Caspase 14

