

TACE

(A disintegrin and metalloproteinase domain 17, TNF-alpha-converting enzyme, TNF-alpha convertase, Snake venom-like protease, [CD156b antigen](#))

- **CAT #:** TACE
- **Enzyme:** TACE, human recombinant (residues 1-671) expressed in *insect cell* with His tag
- **Substrate:** (5-FAM/QXL™) FRET peptide
- **References Compound IC50s:**

<u>Compounds</u>	<u>NNGH</u>	<u>Actinonin</u>	<u>GM6001</u>	<u>MMP Inhibitor III</u>	<u>MMP-9/MMP-13 Inhibitor I</u>
IC50(M)	6.88E-07	1.14E-06	3.03E-08	2.22E-08	1.99E-09

- **Scientific Information:** [P78536](#)
- FAM = Carboxyfluorescein

Control Inhibitor IC50s for TACE

