

---

# TYRO3

Product code: 08-109

Human TYRO3, cytoplasmic domain of [453-890(end) amino acids of accession number NP\_006284.2] was expressed as N-terminal GST fusion protein (76 kDa) using baculovirus expression system.

Assay platform : MSA  
Substrate : CSKtide  
  
ATP ( $\mu$ M)  $K_m$  app / Bin : 80 / 75  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 1.3  
IC50 at 1 mM ATP (nM) : 2.9

---