

---

# RON(MST1R)

Product code 08-152

Human RON, cytoplasmic domain [979-1400(end) amino acids of accession number NP\_002438.1] was expressed as N-terminal GST-fusion protein (75kDa) using baculovirus expression system. GST-RON was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay  
Substrate : Src tide

ATP ( $\mu$ M)  $K_m$  app /  $B_{in}$  : 27 / 25  
Metal : Mg  
Reference compound : Staurosporine  
IC<sub>50</sub> at ATP  $B_{in}$  (nM) : 37  
IC<sub>50</sub> at 1 mM ATP (nM) : 550

---