

---

---

# MAPKAPK5

Product code: 02-144

Full-length human MAPKAPK5 [1-471(end) amino acids of accession number NP\_003659.2] was co-expressed as N-terminal GST-fusion protein (81 kDa) with human His-tagged p38 $\beta$  [1-364(end) amino acids of accession number NP\_002742.3] and human His-tagged MAP2K6 [1-334(end) amino acids of accession number NP\_002749] using baculovirus expression system.

Assay platform : MSA  
Substrate : GS peptide  
  
ATP ( $\mu$ M) Km app / Bin : 12 / 10  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 320  
IC50 at 1 mM ATP (nM) : n.a.

---

---