

---

---

# MAP3K4

Product code: 07-106

Human MAP3K4, catalytic domain [1312-1608(end) amino acids of accession number NP\_005913.2] was expressed as N-terminal GST-fusion protein (61 kDa) using baculovirus expression system.

Assay platform : MSA  
Substrate : MAP2K6  
Cascade Assay \*  
ATP ( $\mu$ M) : 1000  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at 1 mM ATP (nM) : 100

\* MAP2K6/p38 $\alpha$ /Modified Erktide

---

---