

---

---

## MLK3(MAP3K11)

Product code: 09-017

Human MLK3, catalytic domain [99-398 amino acids of accession number NP\_002410.1] was expressed as N-terminal His-tagged protein (37kDa) using baculovirus expression system.

Assay platform : MSA  
Substrate : MAP2K1  
Cascade Assay \*  
ATP ( $\mu$ M) : 1000  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at 1 mM ATP (nM) : 4.8

\* MAP2K1/Erk2/Modified Erktide

---

---