

---

---

## ACK(TNK2)

Product code: 08-196

Human ACK, catalytic domain [110-476 amino acids of accession number NP\_005772.3] was expressed as N-terminal GST-fusion protein (69 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: WASP peptide
ATP ( $\mu$ M) Km app / Bin	: 97 / 100
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 3.2
IC50 at 1 mM ATP (nM)	: 3.8

---

---