

---

# AKT2

Product code: 01-102

Human AKT2, catalytic domain [120-481(end) amino acids of accession number NP\_001617.1] was co-expressed as N-terminal GST-fusion protein (69 kDa) with His-tagged PDK1 [1-556(end) amino acids of accession number NP\_002604.1] using baculovirus expression system. GST-AKT2 was purified by using glutathione sepharose chromatography.

Assay platform : MSA  
Substrate : Crosstide  
  
ATP ( $\mu$ M) Km app / Bin : 110 / 100  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 5.2  
IC50 at 1 mM ATP (nM) : n.a.

---