

---

# AKT3

Product code: 01-103

Human AKT3, catalytic domain [108-479(end) amino acids of accession number NP\_005456.1] was co-expressed as N-terminal GST-fusion protein (70 kDa) with His-tagged PDK1 [1-556(end) amino acids of accession number NP\_002604.1] using baculovirus expression system. GST-AKT3 was purified by using glutathione sepharose chromatography.

Assay platform : MSA  
Substrate : Crosstide  
  
ATP ( $\mu$ M) Km app / Bin : 54 / 50  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 3.2  
IC50 at 1 mM ATP (nM) : n.a.

---