

---

# AKT1

Product code 01-101

Human AKT1, catalytic domain [104-480(end) amino acids of accession number NP\_005154.1] 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-AKT1 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Crosstide

ATP ( $\mu$ M) Km app / Bin : 31 / 50

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 2.7

IC50 at 1 mM ATP (nM) : 22

---