
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 (μ 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.
