
AKT1

Product code: 01-101

Human AKT1, catalytic domain [104-480(end) amino acids and G478S of accession number NP_005154.2] 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.

Assay platform : MSA
Substrate : Crosstide
ATP (μ 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
