
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 (μ 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.
