
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.

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.

