
SGK

Product code: 01-158

Truncated human SGK [61-431(end) amino acids and S422D of accession number NP_005618.2] was co-expressed as N-terminal GST-fusion protein (68 kDa) with His-tagged PDK1 [1-556(end) amino acids of accession number NP_002604.1] using baculovirus expression system.

Assay platform : MSA
Substrate : SGKtide

ATP (μ M) Km app / Bin : 52 / 50
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 10
IC50 at 1 mM ATP (nM) : 99
