
SGK2

Product code: 01-159

Full-length human SGK2 [1-367(end) amino acids and S356D of accession number NP_733794.1] 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	: 58 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 30
IC50 at 1 mM ATP (nM)	: n.a.

