
SLK

Product code: 07-129

Full-length human SLK [1-1152(end) amino acids and S5N of accession number NP_055535.1] was expressed as N-terminal GST-fusion protein (160 kDa) using baculovirus expression system.

Assay platform	:	MSA
Substrate	:	Moesin-derived peptide
ATP (μ M) Km app / Bin	:	36 / 50
Metal	:	Mg
Reference compound	:	Staurosporine
IC50 at ATP Bin (nM)	:	0.32
IC50 at 1 mM ATP (nM)	:	n.a.

