
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.
