
LOK(STK10)

Product code: 07-315

Full-length human LOK [1-968(end) amino acids of accession number BAA35073.1] was expressed as N-terminal GST-fusion protein using baculovirus expression system. GST-LOK was purified by using glutathione sepharose chromatography. GST-LOK was cleaved by PreScission protease and GST-free LOK (114 kDa) was collected as flow-through fraction from glutathione sepharose chromatography.

Assay platform	: MSA
Substrate	: Moesin-derived peptide
ATP (μ M) Km app / Bin	: 100 / 100
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.49
IC50 at 1 mM ATP (nM)	: n.a.
