
QIK(SNF1LK2)

Product code: 02-129

Full-length human QIK(SNF1LK2) [1-926(end) amino acids of accession number NP_056006.1] was expressed as N-terminal GST-fusion protein (132 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: AMARA peptide
ATP (μ M) Km app / Bin	: 42 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.4
IC50 at 1 mM ATP (nM)	: 2.9