
RSK4(RPS6KA6)

Product code: 01-152

Full-length human RSK4 [1-745(end) amino acids of accession number NP_055311.1] was expressed as N-terminal GST-fusion protein (111 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: S6K peptide(N-FL)
ATP (μ M) Km app / Bin	: 20 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.051
IC50 at 1 mM ATP (nM)	: 0.56
