
RSK3(RPS6KA2)

Product code: 01-151

Full-length human RSK3 [1-733(end) amino acids of accession number NP_066958.2] 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	: 9.9 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.11
IC50 at 1 mM ATP (nM)	: 1.7
