
RSK2(RPS6KA3)

Product code: 01-150

Full-length human RSK2 [1-740(end) amino acids of accession number NP_004577.1] 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	: 14 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.10
IC50 at 1 mM ATP (nM)	: 2.1
