
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. GST-RSK4 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

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
