
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. GST-RSK3 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : S6K peptide(N-FL)

ATP (μ M) K_{mapp} / B_{in} : 9.9 / 10

Metal : Mg

Reference compound : Staurosporine

IC₅₀ at ATP B_{in} (nM) : 0.11

IC₅₀ at 1 mM ATP (nM) : 1.7
