
p70S6K(RPS6KB1)

Product code: 01-154

Human p70S6K, catalytic domain [1-421 amino acids and T412E of accession number NP_003152.1] was expressed as N-terminal GST-fusion protein (75 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: S6K2 peptide
ATP (μ M) Km app / Bin	: 14 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.1
IC50 at 1 mM ATP (nM)	: 9.8

