
p70S6K β (RPS6KB2)

Product code: 01-155

Full-length human p70S6K β [1-482(end) amino acids of accession number NP_003943.2] was expressed as N-terminal GST-fusion protein (81 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: S6K2 peptide
ATP (μ M) Km app / Bin	: 3.3 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.7
IC50 at 1 mM ATP (nM)	: n.a.
