

---

---

# RSK1(RPS6KA1)

Product code: 01-149

Full-length human RSK1 [1-735(end) amino acids of accession number NP\_002944.2] was expressed as N-terminal GST-fusion protein (110 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: S6K peptide(N-FL)
ATP ( $\mu$ M) Km app / Bin	: 21 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.14
IC50 at 1 mM ATP (nM)	: 2.5

---

---