
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 (μ 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 |
