

---

## ***p70S6K $\beta$ (RPS6KB2)***

Product code 01-155

Full-length human p70S6K $\beta$  [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. GST-p70S6K $\beta$  was purified by using glutathione sepharose chromatography

Assay platform : Mobility Shift Assay

Substrate : S6K2 peptide

ATP ( $\mu$ M)  $K_{mapp}$  /  $B_{in}$  : 3.3 / 5

Metal : Mg

Reference compound : Staurosporine

IC<sub>50</sub> at ATP  $B_{in}$  (nM) : 2.7

IC<sub>50</sub> at 1 mM ATP (nM) : n.a.

---