

---

## ***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. GST-p70S6K was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay  
Substrate : S6K2 peptide

ATP ( $\mu$ M)  $K_{mapp}$  /  $B_{in}$  : 14 / 10  
Metal : Mg  
Reference compound : Staurosporine  
IC<sub>50</sub> at ATP  $B_{in}$  (nM) : 2.1  
IC<sub>50</sub> at 1 mM ATP (nM) : 9.8

---