



# RSK2,RPS6KA3, MAPKAP-K1 $\beta$ validated with HTRF® KinEASE

## TECHNICAL NOTE

**Enzyme name** : RSK2,RPS6KA3, MAPKAP-K1 beta

**Substrate** : STK S1

**Buffer additives** : 5 mM MgCl<sub>2</sub>, 1 mM DTT

**[RSK2,RPS6KA3, MAPKAP-K1  $\beta$ ]** : 0.33 ng/ $\mu$ L

**Incubation time at RT** : 30 min

**ATP Km** : 26.2  $\mu$ M

**Staurosporine IC50** : 7.282 nM

