



# RSK1,RPS6KA1, MAPKAP-K1 $\alpha$ validated with HTRF® KinEASE

## TECHNICAL NOTE

**Enzyme name** : RSK1,RPS6KA1, MAPKAP-K1 alpha

**Substrate** : STK S1

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

**[RSK1,RPS6KA1, MAPKAP-K1  $\alpha$ ]** : 1 ng/ $\mu$ L

**Incubation time at RT** : 30 min

**ATP Km** : 22.36  $\mu$ M

**Staurosporine IC50** : 51.95 nM

