



# RSK3,RPS6KA2 validated with HTRF® KinEASE

## TECHNICAL NOTE

**Enzyme name** : RSK3,RPS6KA2

**Substrate** : STK S1

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

**[RSK3,RPS6KA2]** : 0.33 ng/μL

**Incubation time at RT** : 50 min

**ATP Km** : 22.03 μM

**Staurosporine IC<sub>50</sub>** : 6.657 nM

