



# ROCK1

## validated with HTRF® KinEASE

### TECHNICAL NOTE

**Enzyme name :** ROCK1

**Substrate :** STK S2

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

**[ROCK1] :** 0.4 ng/μL

**Incubation time at RT :** 20 min

**ATP Km :** 1.29 μM

**Staurosporine IC<sub>50</sub> :** 29.65 nM

