



# PKAC $\alpha$ ,PRKACA validated with HTRF® KinEASE

## TECHNICAL NOTE

**Enzyme name** : PKAC alpha,PRKACA

**Substrate** : STK S2

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

**[PKAC  $\alpha$ ,PRKACA]** : 0.00026 ng/ $\mu$ L

**Incubation time at RT** : 30 min

**ATP Km** : 1.7  $\mu$ M

**Staurosporine IC<sub>50</sub>** : 10.8 nM

