



# PKC $\gamma$ , PRKCG validated with HTRF® KinEASE

## TECHNICAL NOTE

**Enzyme name** : PKC gamma, PRKCG

**Substrate** : STK S1

**Buffer additives** : 10 mM MgCl<sub>2</sub>, 1 mM DTT, lipid activator 1X

**[PKC  $\gamma$ , PRKCG]** : 1 ng/ $\mu$ L

**Incubation time at RT** : 30 min

**ATP Km** : 4.333  $\mu$ M

**Staurosporine IC50** : 6.157 nM

