



# PKC $\zeta$ ,PRKCZ

## validated with HTRF® KinEASE

### TECHNICAL NOTE

**Enzyme name** : PKC zeta,PRKCZ

**Substrate** : STK S1

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

**[PKC  $\zeta$ ,PRKCZ]** : 0.333 ng/ $\mu$ L

**Incubation time at RT** : 30 min

**ATP Km** : 3.254  $\mu$ M

**Staurosporine IC50** : 119 nM

