



# CGK2,PRKG2 validated with HTRF® KinEASE

## TECHNICAL NOTE

**Enzyme name :** CGK2,PRKG2  
**Substrate :** STK S2  
**Buffer additives :** 5 mM MgCl<sub>2</sub>, 1 mM DTT  
**[CGK2,PRKG2] :** 0.037 ng/μL  
**Incubation time at RT :** 40 min  
**ATP Km :** 79.32 μM  
**Staurosporine IC<sub>50</sub> :** 5.52 nM

