



# PLK4

## validated with HTRF® KinEASE

### TECHNICAL NOTE

**Enzyme name** : PLK4

**Substrate** : STK S2

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

**[PLK4]** : 1 ng/μL

**Incubation time at RT** : 60 min

**ATP Km** : 26 μM

**Staurosporine IC<sub>50</sub>** : 75 nM

