



# smMLCK,MYLK validated with HTRF® KinEASE

## TECHNICAL NOTE

**Enzyme name** : smMLCK,MYLK

**Substrate** : STK S1

**Buffer additives** : 10 mM MgCl<sub>2</sub>, 1 mM DTT, 500 μM CaCl<sub>2</sub>, 1 μM Calmodulin

**[smMLCK,MYLK]** : 1 ng/μL

**Incubation time at RT** : 50 min

**ATP Km** : n.d. μM

**Staurosporine IC<sub>50</sub>** : 17.25 nM

