



# DCAMKL2

## validated with HTRF® KinEASE

### TECHNICAL NOTE

**Enzyme name :** DCAMKL2

**Substrate :** STK S1

**Buffer additives :** 5 mM MgCl<sub>2</sub>, 1 mM DTT, 500 μM CaCl<sub>2</sub>, 1 μM Calmodulin

**[DCAMKL2] :** 0.037 ng/μL

**Incubation time at RT :** 50 min

**ATP Km :** 22.82 μM

**Staurosporine IC<sub>50</sub> :** 90.1 nM

