



# EPHA1

## validated with HTRF® KinEASE

### TECHNICAL NOTE

**Enzyme name :** EPHA1

**Substrate :** TK

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

**[EPHA1] :** 0.037 ng/μL

**Incubation time at RT :** 50 min

**ATP Km :** 1 μM

**Staurosporine IC50 :** 1486 nM

