



# EPHA2

## validated with HTRF® KinEASE

### TECHNICAL NOTE

**Enzyme name :** EPHA2

**Substrate :** TK

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

**[EPHA2] :** 0.037 ng/μL

**Incubation time at RT :** 50 min

**ATP Km :** 6.3 μM

**Staurosporine IC50 :** 4495 nM

