



# EPHA4

## validated with HTRF® KinEASE

### TECHNICAL NOTE

**Enzyme name :** EPHA4

**Substrate :** TK

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

**[EPHA4] :** 0.037 ng/μL

**Incubation time at RT :** 50 min

**ATP Km :** 5.9 μM

**Staurosporine IC50 :** 5851 nM

