



# EPHB4

## validated with HTRF® KinEASE

### TECHNICAL NOTE

**Enzyme name :** EPHB4

**Substrate :** TK

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

**[EPHB4] :** 0.0123 ng/μL

**Incubation time at RT :** 20 min

**ATP Km :** 0.93 μM

**Staurosporine IC50 :** 1743 nM

