



# EPHB2

## validated with HTRF® KinEASE

### TECHNICAL NOTE

**Enzyme name :** EPHB2

**Substrate :** TK

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

**[EPHB2] :** 0.0123 ng/μL

**Incubation time at RT :** 30 min

**ATP Km :** 1.1 μM

**Staurosporine IC<sub>50</sub> :** 112.1 nM

