



# FGFR4

## validated with HTRF® KinEASE

### TECHNICAL NOTE

**Enzyme name :** FGFR4

**Substrate :** TK

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

**[FGFR4] :** 0.333 ng/μL

**Incubation time at RT :** 40 min

**ATP Km :** 113 μM

**Staurosporine IC<sub>50</sub> :** 251.2 nM

