



# FES,Fps validated with HTRF® KinEASE

## TECHNICAL NOTE

**Enzyme name :** FES,Fps

**Substrate :** TK

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

**[FES,Fps] :** 0.14 ng/μL

**Incubation time at RT :** 50 min

**ATP Km :** 17.9 μM

**Staurosporine IC50 :** 310 nM

