



# FRK,PTK5 validated with HTRF® KinEASE

## TECHNICAL NOTE

**Enzyme name :** FRK,PTK5

**Substrate :** TK

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

**[FRK,PTK5] :** 0.037 ng/μL

**Incubation time at RT :** 50 min

**ATP Km :** 1.887 μM

**Staurosporine IC<sub>50</sub> :** 24.1 nM

