



# FLT3

## validated with HTRF® KinEASE

### TECHNICAL NOTE

**Enzyme name :** FLT3

**Substrate :** TK

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

**[FLT3] :** 1 ng/μL

**Incubation time at RT :** 40 min

**ATP Km :** 4.29 μM

**Staurosporine IC50 :** 29.2 nM

