



# ZAP70

## validated with HTRF® KinEASE

### TECHNICAL NOTE

**Enzyme name :** ZAP70

**Substrate :** TK

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

**[ZAP70] :** 0.037 ng/μL

**Incubation time at RT :** 50 min

**ATP Km :** 0.57 μM

**Staurosporine IC<sub>50</sub> :** 16.93 nM

