



# RON,MST1R validated with HTRF® KinEASE

## TECHNICAL NOTE

**Enzyme name :** RON,MST1R

**Substrate :** TK

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

**[RON,MST1R] :** 0.111 ng/μL

**Incubation time at RT :** 50 min

**ATP Km :** 2.3 μM

**Staurosporine IC<sub>50</sub> :** 2306 nM

