



# PIM1

## validated with HTRF® KinEASE

### TECHNICAL NOTE

**Enzyme name :** PIM1

**Substrate :** STK S3

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

**[PIM1] :** 0.33 ng/μL

**Incubation time at RT :** 50 min

**ATP Km :** 36.35 μM

**Staurosporine IC<sub>50</sub> :** 24.71 nM

