



# PDGFR $\beta$ , PDGFRB validated with HTRF® KinEASE

## TECHNICAL NOTE

**Enzyme name :** PDGFR beta, PDGFRB

**Substrate :** TK

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

**[PDGFR  $\beta$ , PDGFRB] :** 0.333 ng/ $\mu$ L

**Incubation time at RT :** 50 min

**ATP Km :** 27.3  $\mu$ M

**Staurosporine IC50 :** 48.6 nM

