



PDGFR α , PDGFRA validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : PDGFR alpha, PDGFRA

Substrate : TK

Buffer additives : 5 mM MgCl₂, 1 mM MnCl₂, 1 mM DTT, SEB

[PDGFR α , PDGFRA] : 0.037 ng/ μ L

Incubation time at RT : 20 min

ATP Km : 0.49 μ M

Staurosporine IC50 : 2.15 nM

