



FAK,PTK2 validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : FAK,PTK2

Substrate : TK

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

[FAK,PTK2] : 0.111 ng/μL

Incubation time at RT : 50 min

ATP Km : 13.75 μM

Staurosporine IC50 : 119.2 nM

