



PYK2,PTK2B validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : PYK2,PTK2B

Substrate : TK

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

[PYK2,PTK2B] : 0.037 ng/μL

Incubation time at RT : 30 min

ATP Km : 1.1 μM

Staurosporine IC₅₀ : 2.173 nM

