



PRK2,PKN2 validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : PRK2,PKN2

Substrate : STK S1

Buffer additives : 5 mM MgCl₂, 1 mM DTT

[PRK2,PKN2] : 0.33 ng/μL

Incubation time at RT : 50 min

ATP Km : 34.9 μM

Staurosporine IC₅₀ : 20.42 nM

