



DAPK3,ZIPK validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : DAPK3,ZIPK

Substrate : STK S1

Buffer additives : 10 mM MgCl₂, 1 mM DTT

[DAPK3,ZIPK] : 1 ng/μL

Incubation time at RT : 50 min

ATP Km : 1.14 μM

Staurosporine IC₅₀ : 3.2 nM

