



DAPK1

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : DAPK1

Substrate : STK S1

Buffer additives : 10 mM MgCl₂, 1 mM DTT

[DAPK1] : 0.11 ng/μL

Incubation time at RT : 30 min

ATP Km : 0.1375 μM

Staurosporine IC₅₀ : 7.742 nM

