



DDR2

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : DDR2

Substrate : TK

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

[DDR2] : 0.111 ng/μL

Incubation time at RT : 50 min

ATP Km : 2.1 μM

Staurosporine IC50 : 97.21 nM

