



DYRK2

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : DYRK2

Substrate : STK S3

Buffer additives : 5 mM MgCl₂, 1 mM DTT, 2 mg/ml Casein

[DYRK2] : 1 ng/μL

Incubation time at RT : 50 min

ATP Km : 17.41 μM

Staurosporine IC50 : 26.98 nM

