



CaMK1 α ,CAMK1 validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : CaMK1 alpha,CAMK1

Substrate : STK S1

Buffer additives : 5 mM MgCl₂, 1 mM DTT, 500 μ M CaCl₂, 1 μ M Calmodulin

[CaMK1 α ,CAMK1] : 0.0123 ng/ μ L

Incubation time at RT : 50 min

ATP Km : 31.66 μ M

Staurosporine IC50 : 23.89 nM

