



CaMK2 β ,CAMK2B validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : CaMK2 beta,CAMK2B

Substrate : STK S1

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

[CaMK2 β ,CAMK2B] : 0.00137 ng/ μ L

Incubation time at RT : 50 min

ATP Km : 5.398 μ M

Staurosporine IC50 : 3.171 nM

