



CaMK2 δ , CAMK2D validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : CaMK2 delta, CAMK2D

Substrate : STK S1

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

[CaMK2 δ , CAMK2D] : 0.00137 ng/ μ L

Incubation time at RT : 20 min

ATP Km : 0.9511 μ M

Staurosporine IC50 : 2.647 nM

