



CaMK2 γ ,CAMK2G validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : CaMK2 gamma,CAMK2G

Substrate : STK S1

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

[CaMK2 γ ,CAMK2G] : 0.0123 ng/ μ L

Incubation time at RT : 50 min

ATP Km : 1.635 μ M

Staurosporine IC50 : 2.627 nM

