



PKG1 α ,PRKG1 validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : PKG1 alpha,PRKG1

Substrate : STK S2

Buffer additives : 5 mM MgCl₂, 1 mM DTT, 10 μ M cGMP

[PKG1 α ,PRKG1] : 0.0123 ng/ μ L

Incubation time at RT : 30 min

ATP Km : 0.8219 μ M

Staurosporine IC50 : 13 nM

