



AMPK α 2/ β 2/ γ 1,PRKA A2/B2/G1 validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : AMPK alpha2/beta2/gamma1,PRKA A2/B2/G1

Substrate : STK S1

Buffer additives : 5 mM MgCl₂, 1 mM DTT, 50 μ M AMP

[AMPK α 2/ β 2/ γ 1,PRKA A2/B2/G1] : 0.004 ng/ μ L

Incubation time at RT : 50 min

ATP Km : 17.92 μ M

Staurosporine IC50 : 0.18 nM

