



AKT1,PKB α validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : AKT1,PKB alpha

Substrate : STK S3

Buffer additives : 5 mM MgCl₂, 1 mM DTT

[AKT1,PKB α] : 0.0014 ng/ μ L

Incubation time at RT : 50 min

ATP Km : 25.28 μ M

Staurosporine IC₅₀ : 17.18 nM

