



AKT3,PKB γ validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : AKT3,PKB gamma

Substrate : STK S3

Buffer additives : 5 mM MgCl₂, 1 mM DTT

[AKT3,PKB γ] : 0.0014 ng/ μ L

Incubation time at RT : 50 min

ATP Km : 18.39 μ M

Staurosporine IC₅₀ : 29.06 nM

