



AKT2,PKB β validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : AKT2,PKB beta

Substrate : STK S3

Buffer additives : 5 mM MgCl₂, 1 mM DTT

[AKT2,PKB β] : 0.0041 ng/ μ L

Incubation time at RT : 50 min

ATP Km : 88.26 μ M

Staurosporine IC₅₀ : 39.53 nM

