



TSSK1,STK22D validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : TSSK1,STK22D

Substrate : STK S1

Buffer additives : 10 mM MgCl₂, 1 mM DTT

[TSSK1,STK22D] : 0.11 ng/μL

Incubation time at RT : 50 min

ATP Km : 8.36 μM

Staurosporine IC₅₀ : 0.23 nM

