



TRKA,NTRK1 validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : TRKA,NTRK1

Substrate : TK

Buffer additives : 5 mM MgCl₂, 1 mM DTT

[TRKA,NTRK1] : 0.111 ng/μL

Incubation time at RT : 30 min

ATP Km : 14.68 μM

Staurosporine IC₅₀ : 7.45 nM

