



TRKC, NTRK3 validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : TRKC, NTRK3

Substrate : TK

Buffer additives : 5 mM MgCl₂, 1 mM DTT

[TRKC, NTRK3] : 0.037 ng/μL

Incubation time at RT : 40 min

ATP Km : 25.64 μM

Staurosporine IC₅₀ : 315 nM

