



# TRKB,NTRK2 validated with HTRF® KinEASE

## TECHNICAL NOTE

**Enzyme name :** TRKB,NTRK2

**Substrate :** TK

**Buffer additives :** 5 mM MgCl<sub>2</sub>, 1 mM MnCl<sub>2</sub>, 1 mM DTT

**[TRKB,NTRK2] :** 0.037 ng/μL

**Incubation time at RT :** 40 min

**ATP Km :** 4.77 μM

**Staurosporine IC<sub>50</sub> :** 5.15 nM

