



BRK,PTK6 validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : BRK,PTK6

Substrate : TK

Buffer additives : 5 mM MgCl₂, 1 mM MnCl₂, 1 mM DTT, SEB

[BRK,PTK6] : 0.333 ng/μL

Incubation time at RT : 60 min

ATP Km : 31.9 μM

Staurosporine IC₅₀ : 38050 nM

