



EGFR

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : EGFR

Substrate : TK

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

[EGFR] : 0.0041 ng/μL

Incubation time at RT : 40 min

ATP Km : 1.57 μM

Staurosporine IC₅₀ : 242 nM

