



TYK2

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : TYK2

Substrate : TK

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

[TYK2] : 1 ng/μL

Incubation time at RT : 50 min

ATP Km : 5.06 / 10 μM

Staurosporine IC₅₀ : 117.6 nM

