



CSK

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : CSK

Substrate : TK

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

[CSK] : 0.333 ng/μL

Incubation time at RT : 40 min

ATP Km : 4.39 μM

Staurosporine IC₅₀ : 468 nM

