



FMS,CSF1R validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : FMS,CSF1R

Substrate : TK

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

[FMS,CSF1R] : 0.0123 ng/μL

Incubation time at RT : 40 min

ATP Km : 2.4 μM

Staurosporine IC50 : 4.978 nM

