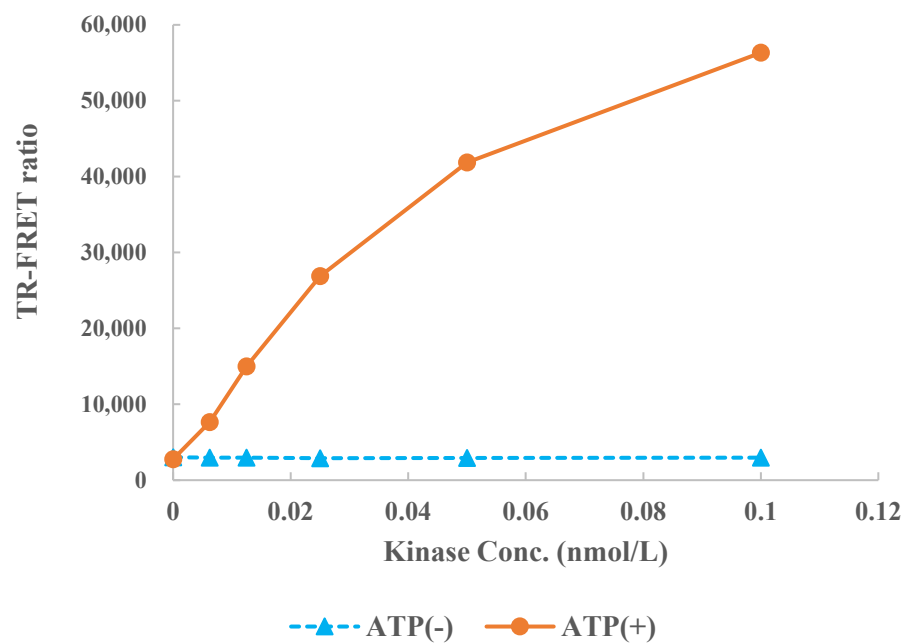


ALK4(ACVR1B)



Kinase	ALK4(ACVR1B), 09-135
Substrate	ULight-TopoII α peptide
Substrate Manufacturer	Revvity, TRF0130-D
Substrate Conc.	50 nmol/L
Antibody	LANCE Eu-anti-Topo II- α (Thr1342) antibody
Antibody Manufacturer	Revvity, TRF0218-D
Antibody Conc.	0.4 nmol/L
ATP Conc.	100 μ mol/L
Metal	Mn
Metal Conc.	5 mmol/L
Reaction Time	1 hour