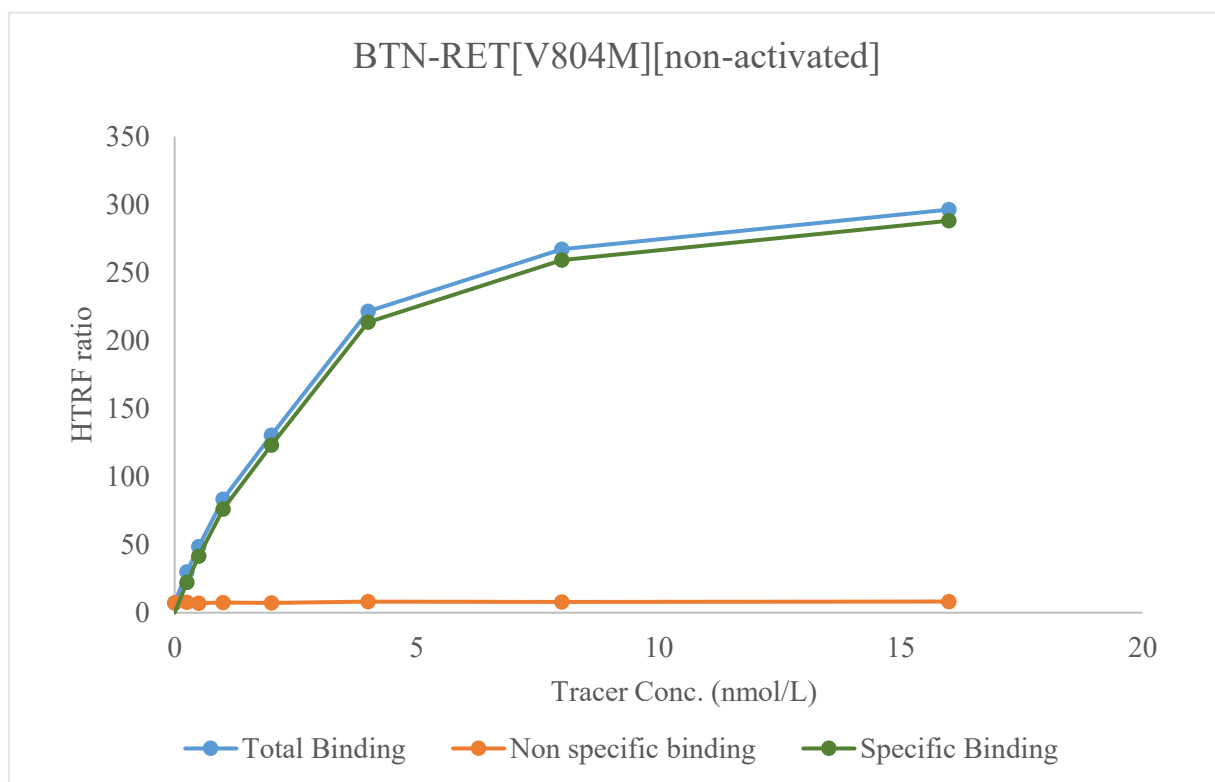


Kinase **BTN-RET[V804M][non-activated]**

Cat.# 08-460-23N



Tracer	Thermo Scientific
	Kinase Tracer 236, PV5592
KD	1.4 nmol/L
Kinase Conc.	5 nmol/L
Eu labeled-streptavidin	5 nmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 Hour