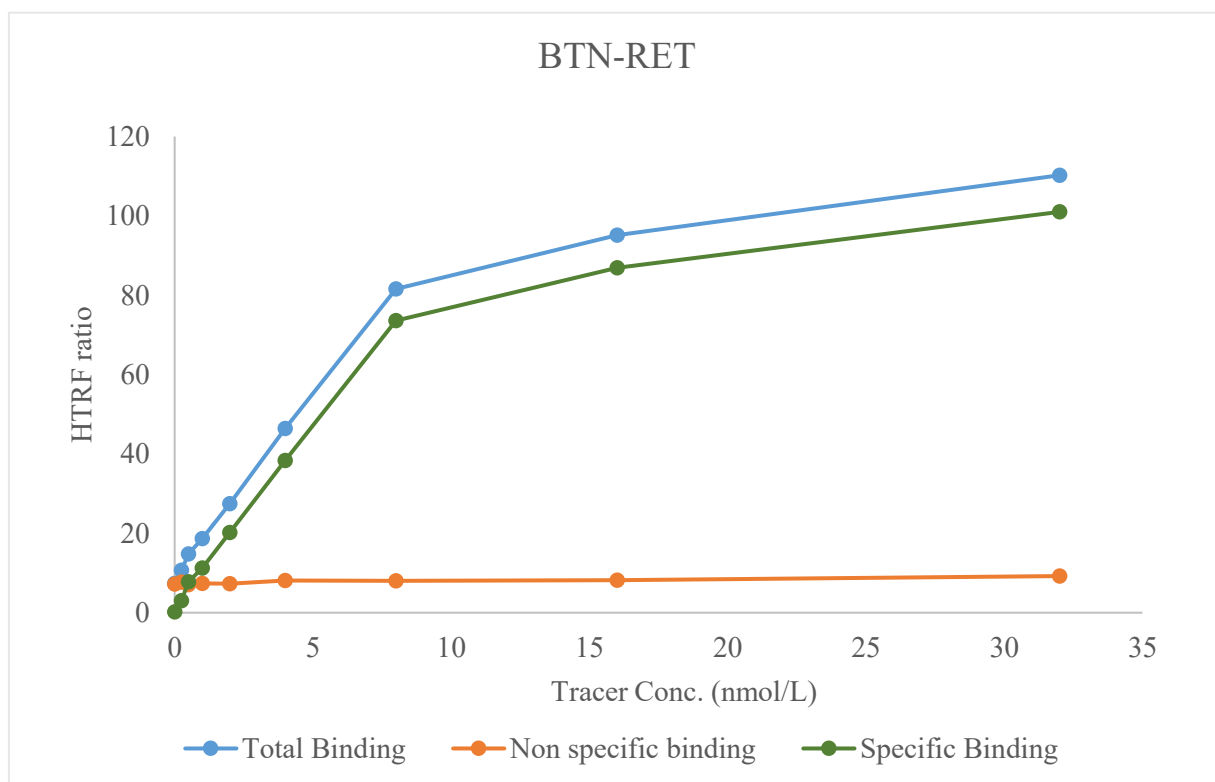


Kinase **BTN-RET**
Cat.# 08-459-20N



Tracer	Thermo Scientific
	Kinase Tracer 236, PV5592
KD	4.6 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