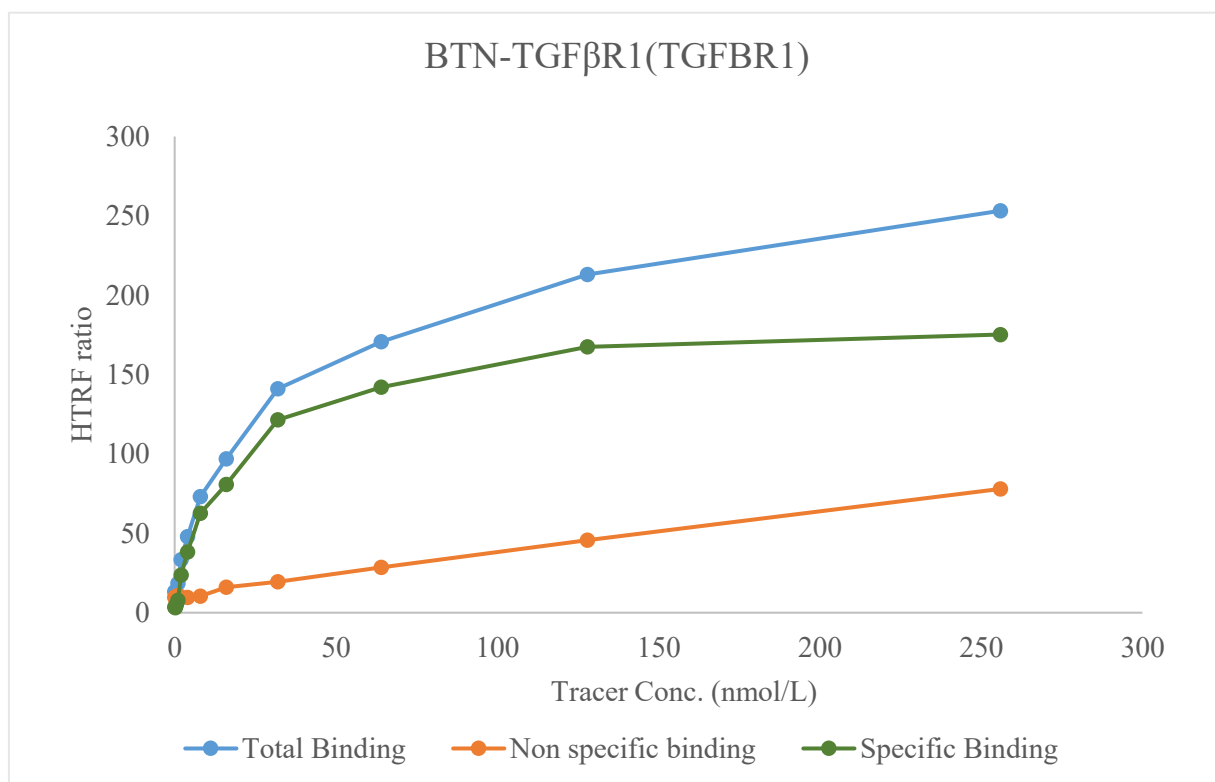


Kinase **BTN-TGF β R1(TGFBR1)**

Cat.# 09-441-20N



Tracer	Thermo Scientific
	Kinase Tracer 178, PV5593
KD	18 nmol/L
Kinase Conc.	5 nmol/L
Eu labeled-streptavidin	5 nmol/L
Metal	Mn
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 Hour