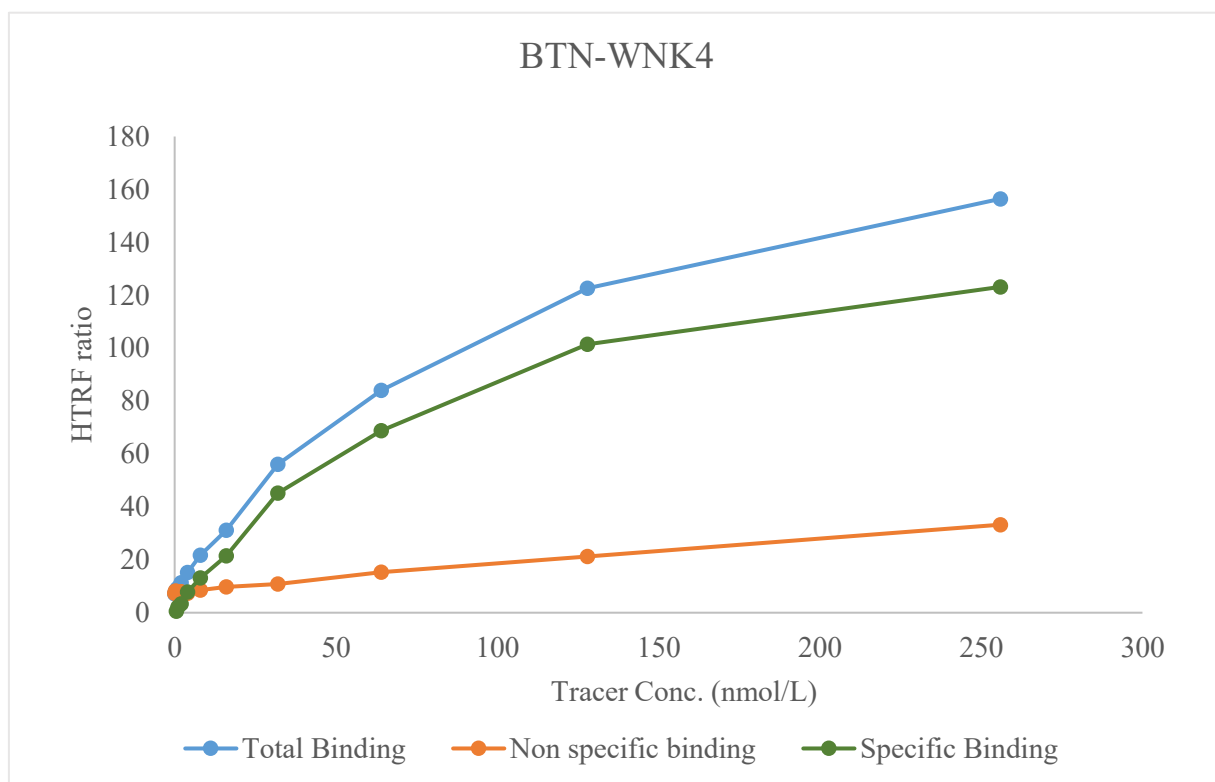


Kinase **BTN-WNK4**

Cat.# 05-482-20N



Tracer	Thermo Scientific
	Kinase Tracer 1710, PV6088
KD	89 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