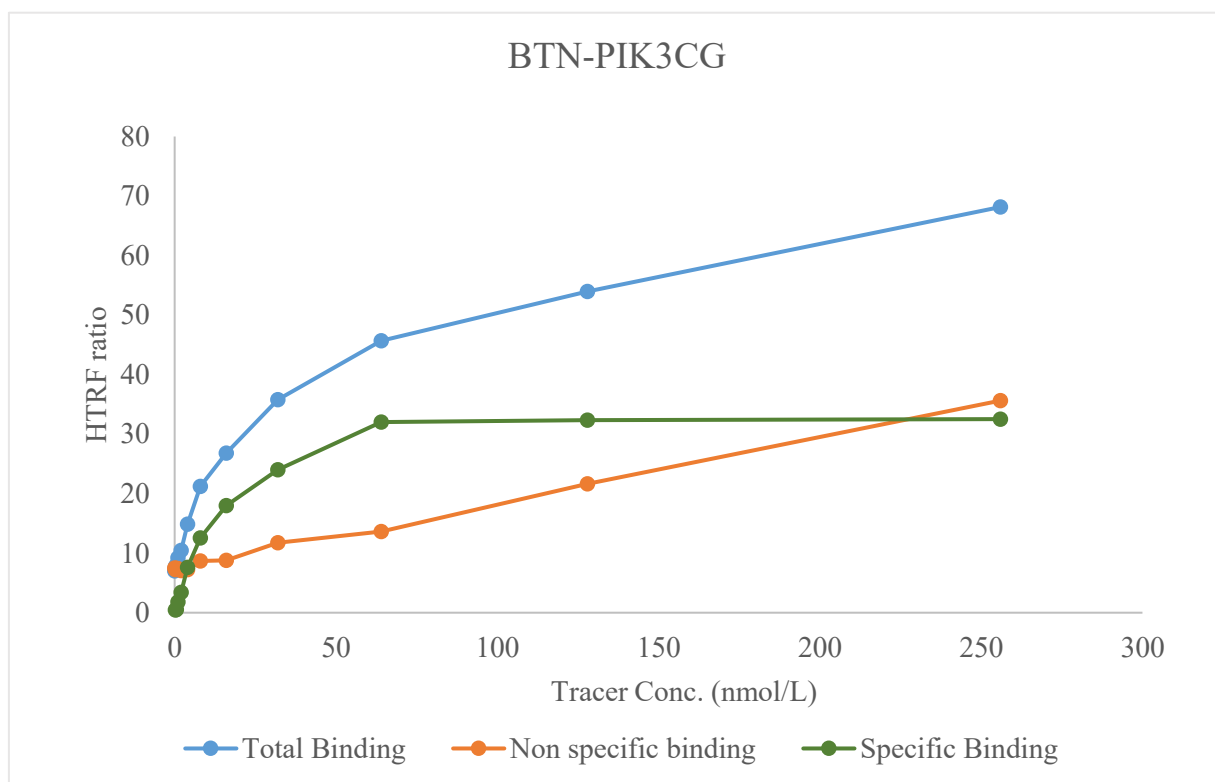


Kinase **BTN-PIK3CG**

Cat.# 11-404-20N



Tracer	Thermo Scientific
	Kinase Tracer 314, PV6087
KD	15 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