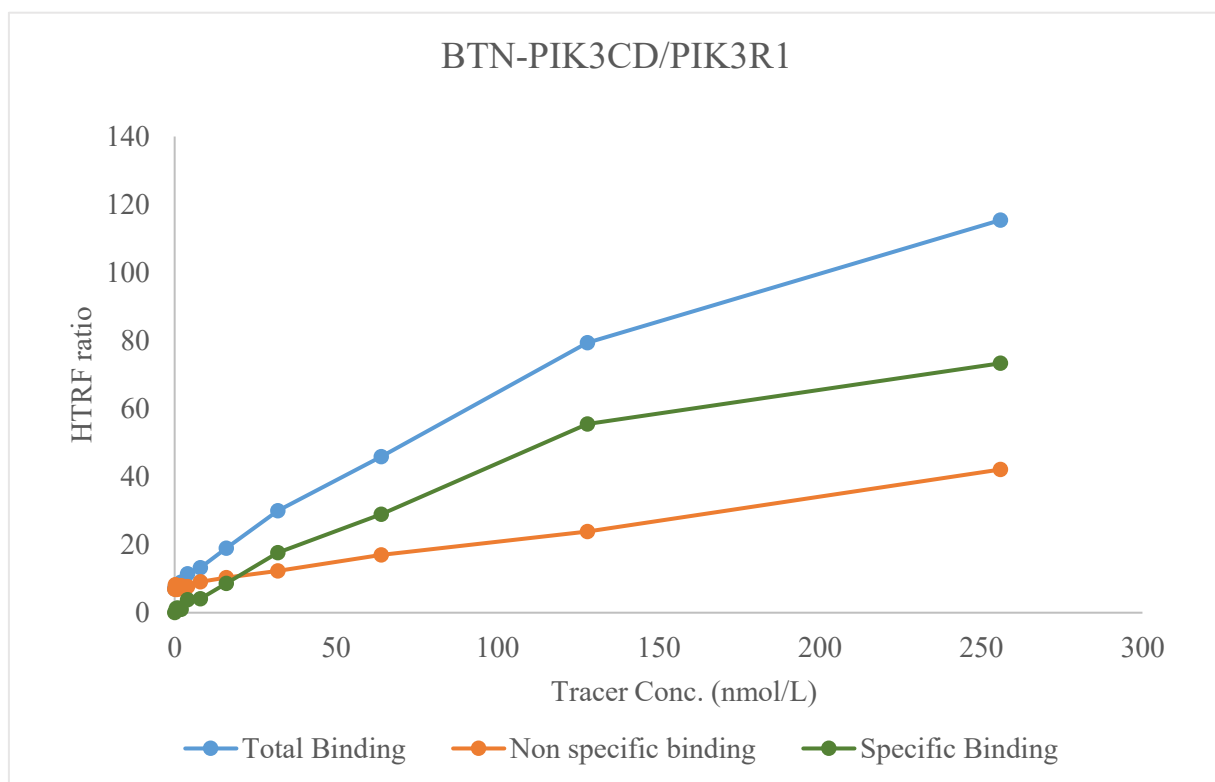


Kinase **BTN-PIK3CD/PIK3R1**

Cat.# 11-403-20N



Tracer	Thermo Scientific
KD	Kinase Tracer 1710, PV6088
	200 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