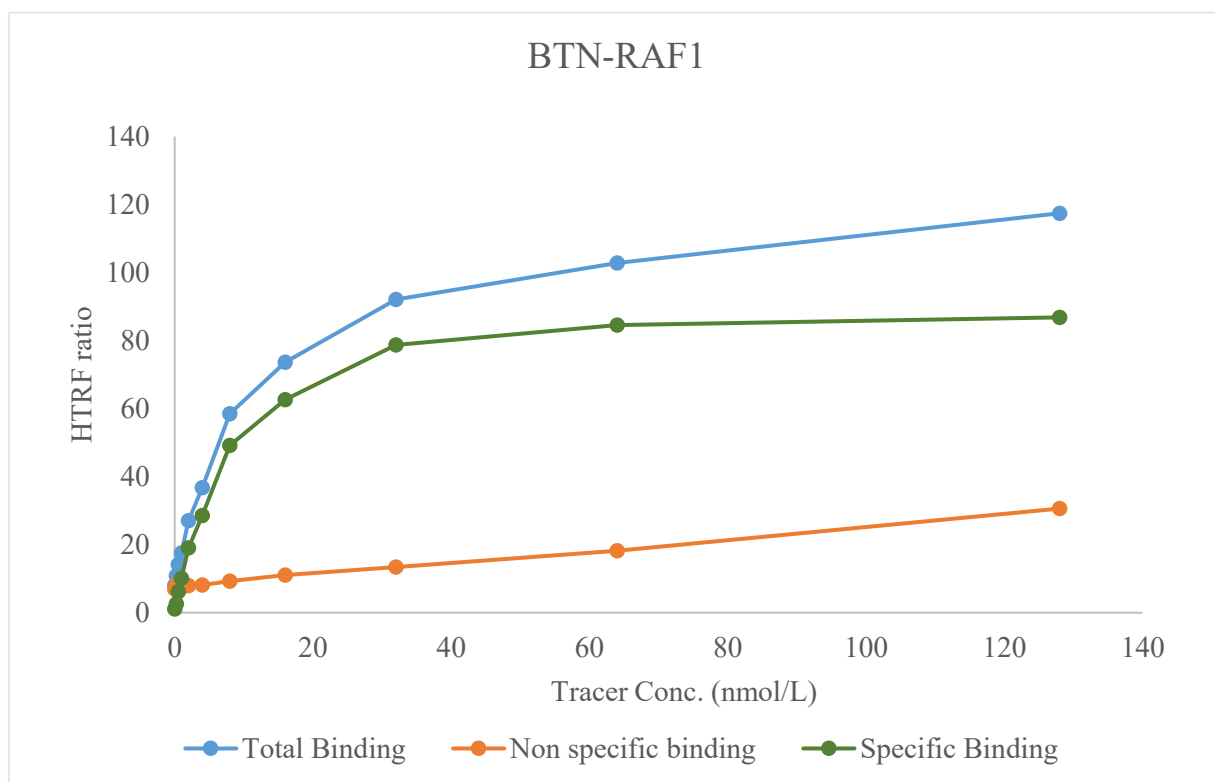


Kinase **BTN-RAF1**

Cat.# 09-425-20N



Tracer	Thermo Scientific
	Kinase Tracer 178, PV5593
KD	8.4 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