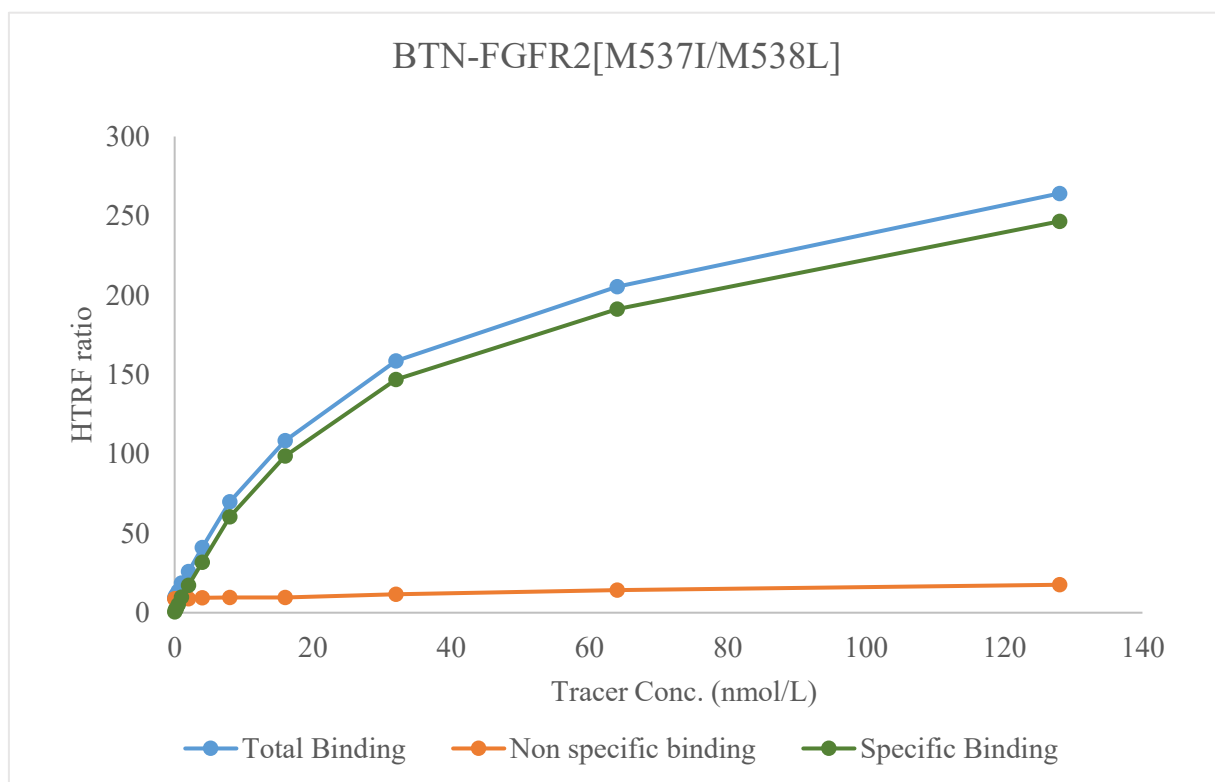


Kinase **BTN-FGFR2[M537I/M538L]**

Cat.# 08-924-20N



Tracer	revvity Inc. Staurosporine-Red, (#62KB01REDC)
KD	35 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