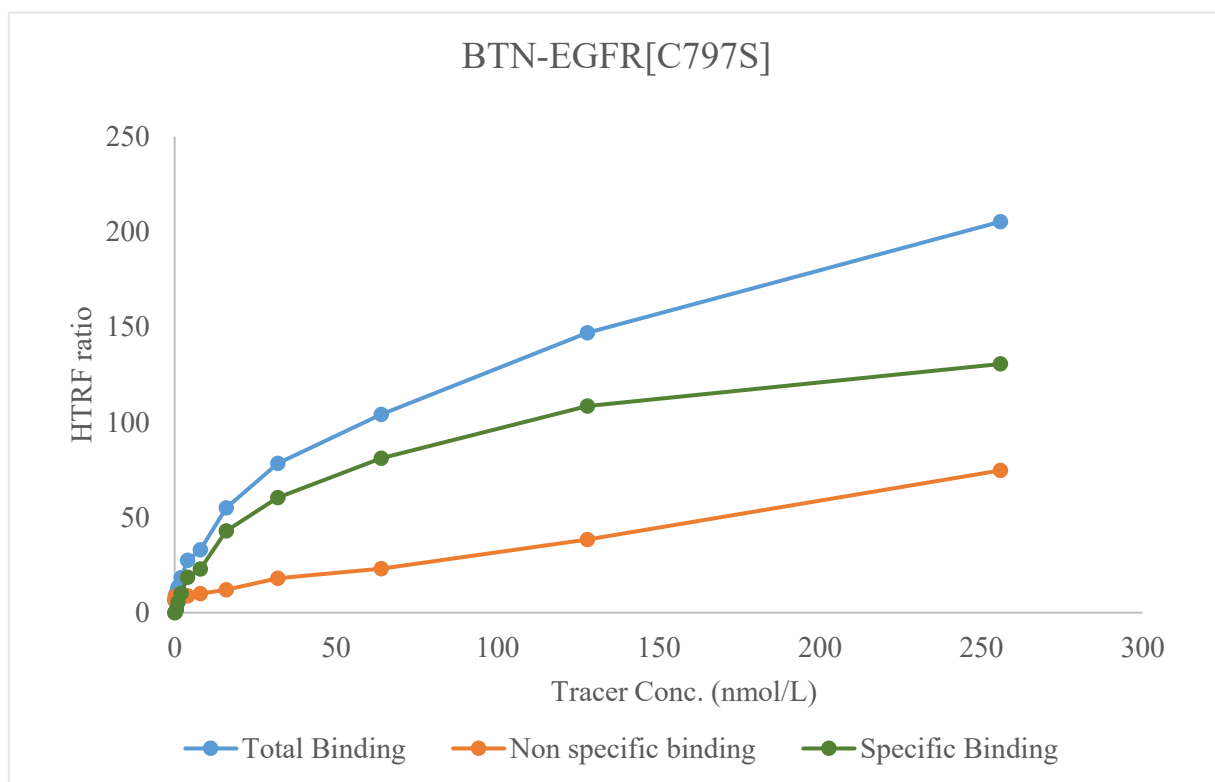


Kinase **BTN-EGFR[C797S]**

Cat.# 08-761-20N



Tracer	revvity Inc. Dasatinib-Red, (#62KB02REDC)
KD	46 nmol/L
Kinase Conc.	5 nmol/L
Eu labeled-streptavidin	5 nmol/L
Metal	Mg, Mn
Metal Conc.	5, 1 mmol/L
Additives	-
Reaction Time	1 Hour