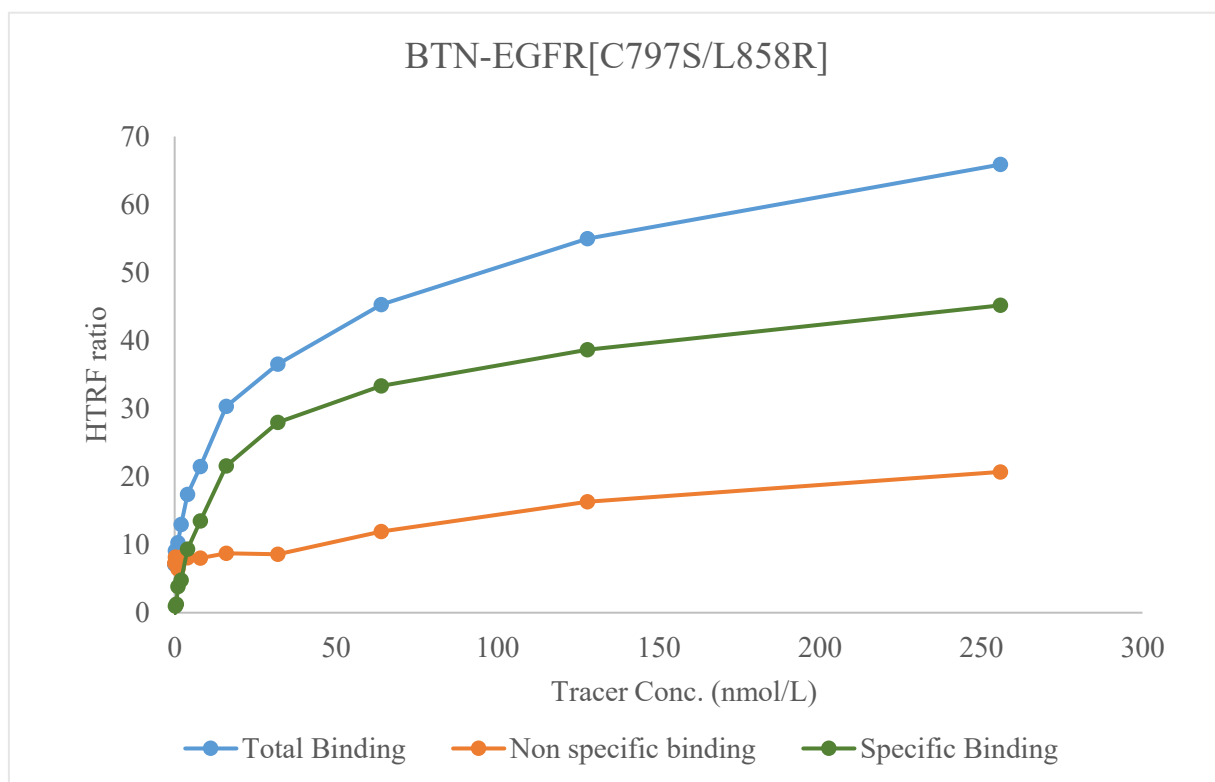


Kinase **BTN-EGFR[C797S/L858R]**

Cat.# 08-763-20N



Tracer	revvity Inc. Staurosporine-Red, (#62KB01REDC)
KD	19 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