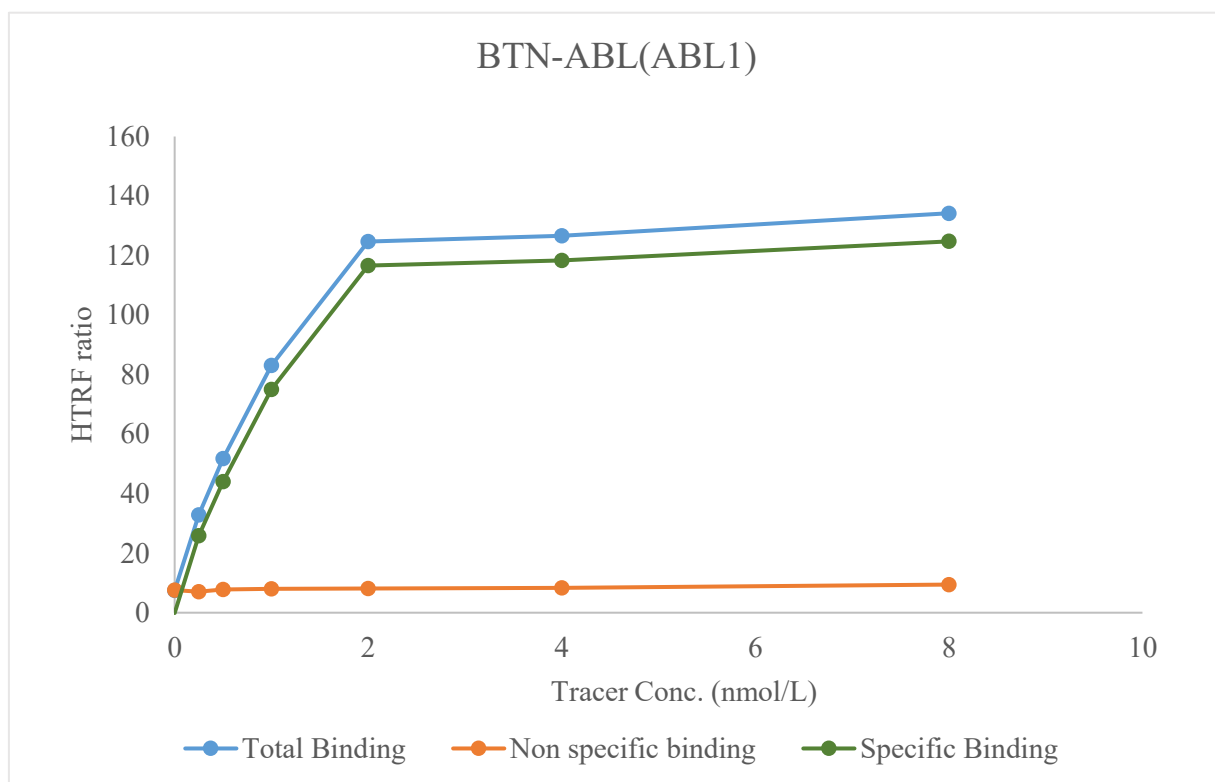


Kinase **BTN-ABL(ABL1)**

Cat.# 08-401-20N



Tracer	revvity Inc. Dasatinib-Red, (#62KB02REDC)
KD	0.76 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