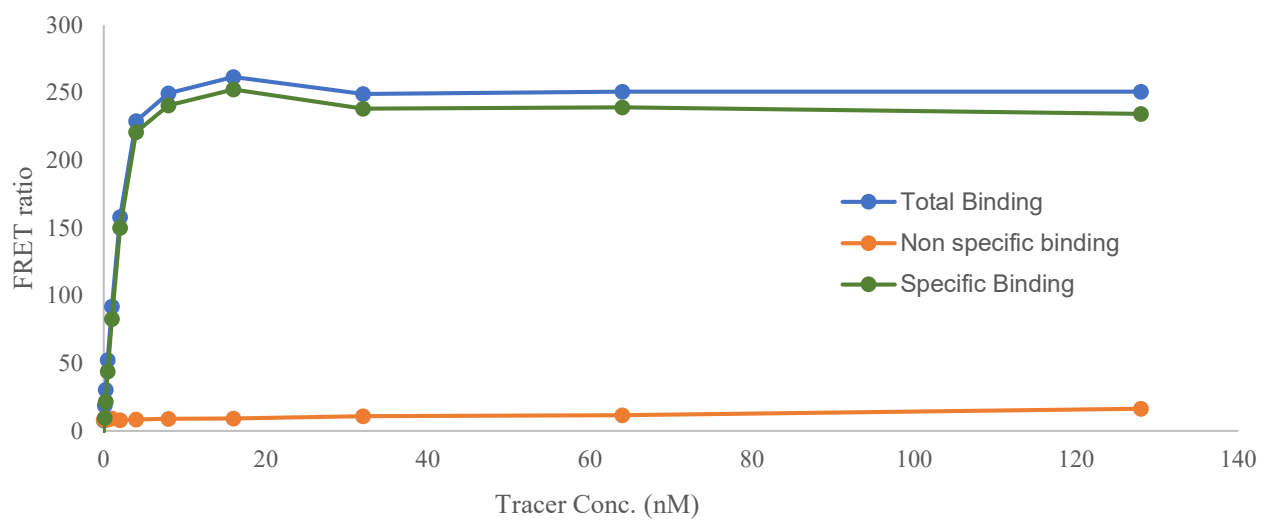


Kinase **BTN-EGFR[d746-750/T790M/C797S/L858R]**
 Cat.# 08-711-20N
 Tracer Revvity Inc.
 Staurosporine-Red (#62KB01REDC)

BTN-EGFR[d746-750/T790M/C797S/L858R]



	KD (nM)	1.61
Kinase conc.		5nM
Eu labeled-streptavidin		5nM

Follow the guide to set up your HTRF® binding assay

Determination of the dissociation constant (Kd) of the fluorescent tracer using HTRF® kinase binding

