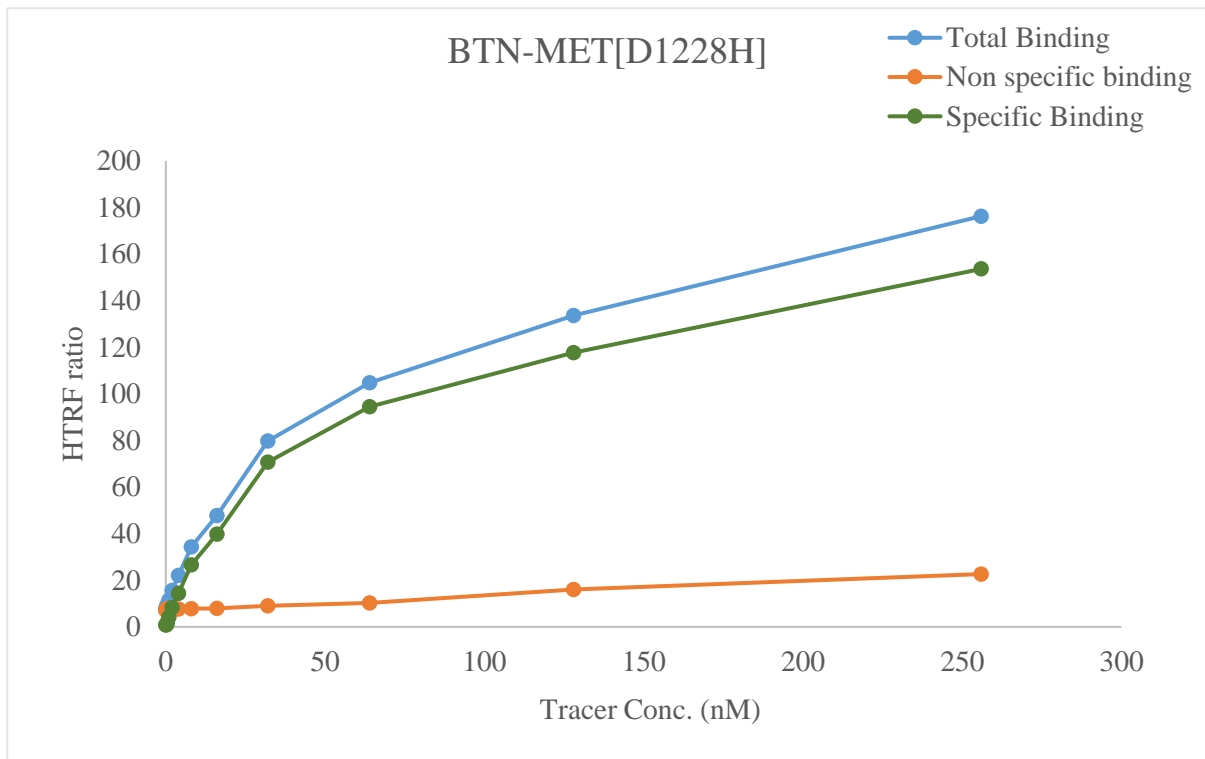


Kinase **BTN-MET[D1228H]**
 Cat.# 08-440-20N
 Tracer revvity Inc.
 Staurosporine-Red (#62KB01REDC)



KD (nM) 55

Kinase conc. 5 nM
 Eu labeled-streptavidin 5 nM

Follow the guide to set up your HTRF[®] binding assay

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

