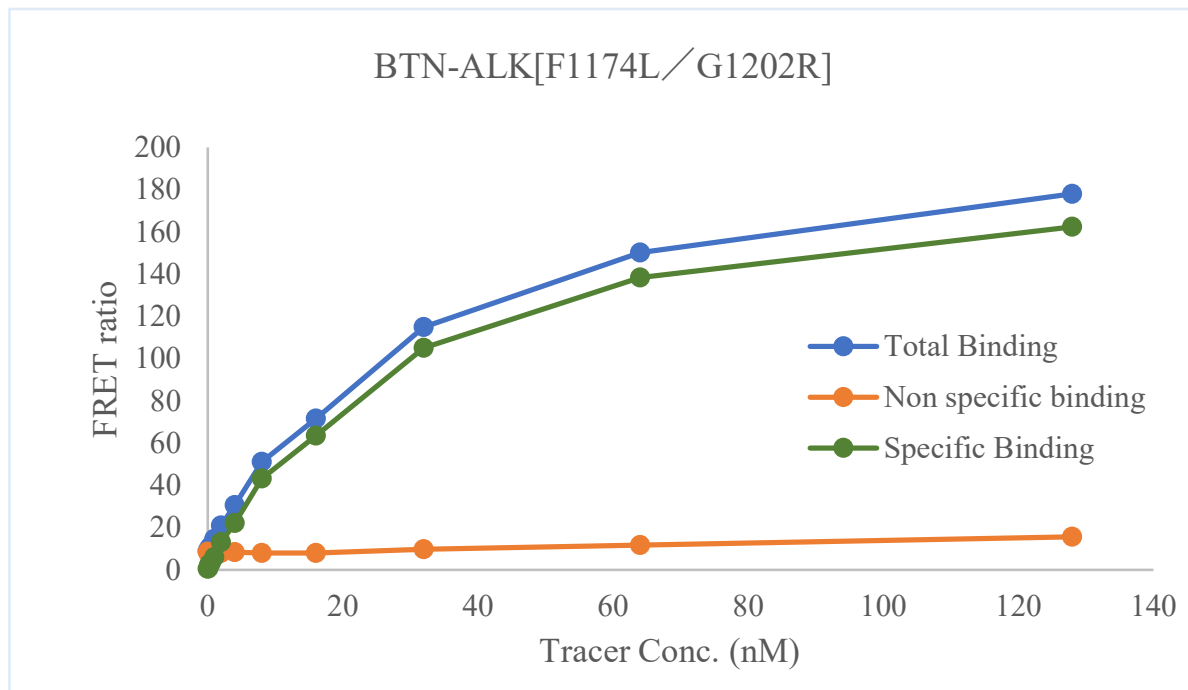


Kinase **BTN-ALK[F1174L/G1202R]**

Cat.# 08-768-20N

Tracer Revvity Inc.

Staurosporine-Red



KD (nM) 32.78

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

