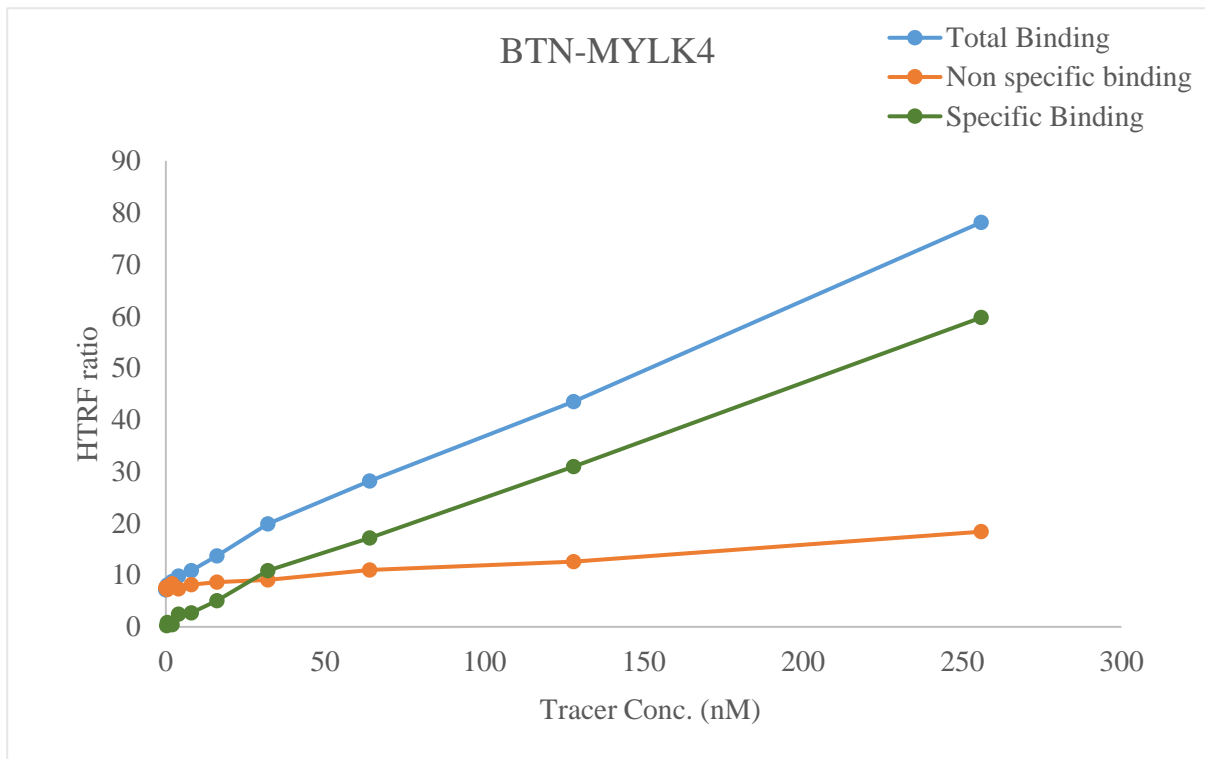


Kinase **BTN-MYLK4**

Cat.# 02-461-20N

Tracer PerkinElmer Inc.

Sunitinib-Red (#62KB03REDC)



KD (nM) ; Not Determined

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

