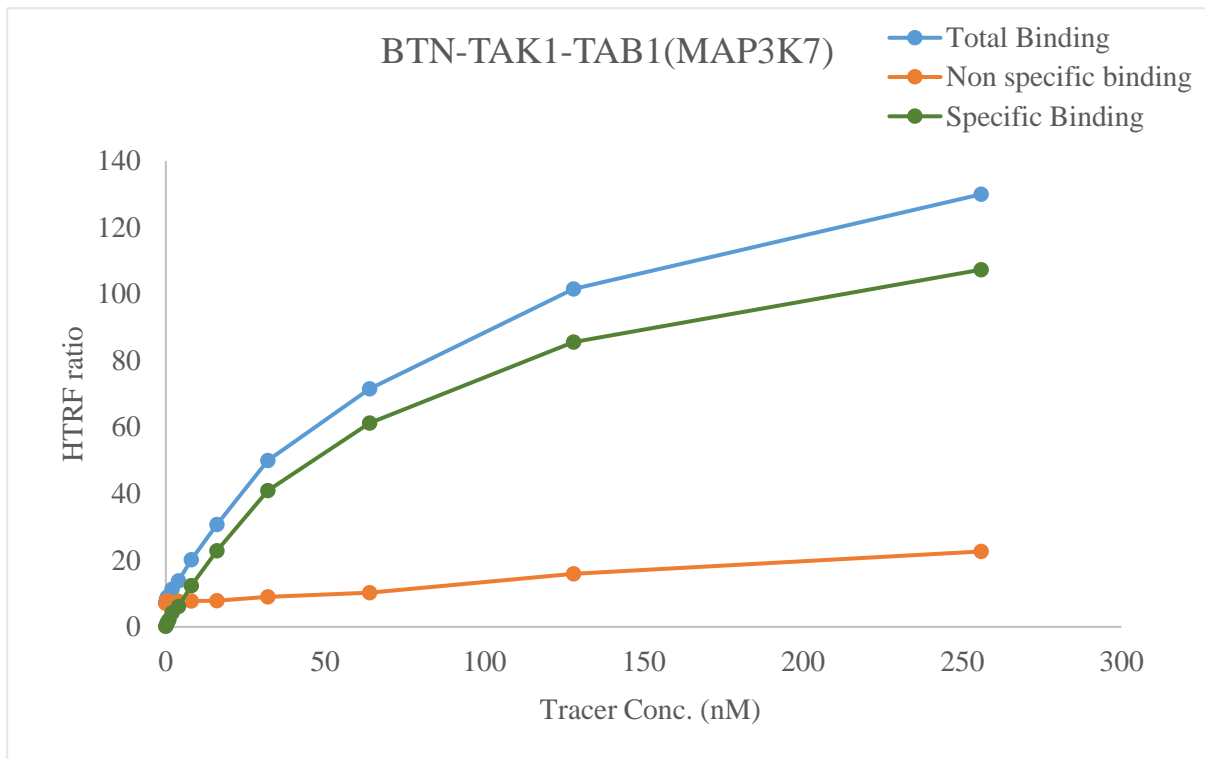


Kinase **BTN-TAK1-TAB1(MAP3K7)**

Cat.# 09-419-21C

Tracer revvity Inc.

Staurosporine-Red (#62KB01REDC)



KD (nM) 84

Kinase conc. 5 nM

Eu labeled-streptavidin 5 nM

Follow the guide to set up your HTRF<sup>®</sup> binding assay

Determination of the dissociation constant (Kd) of the fluorescent tracer using HTRF<sup>®</sup> kinase binding

