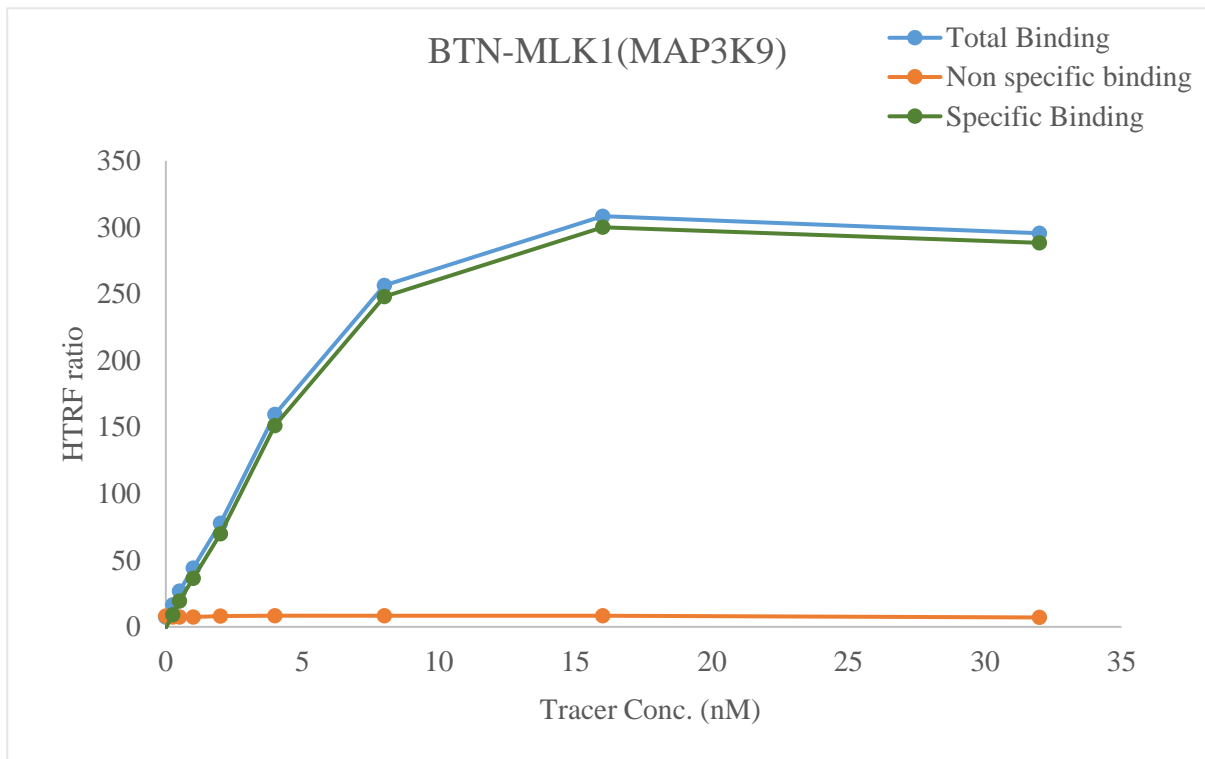


Kinase **BTN-MLK1(MAP3K9)**

Cat.# 09-415-20N

Tracer PerkinElmer Inc.

Staurosporine-Red (#62KB01REDC)



KD (nM) 2.2

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

