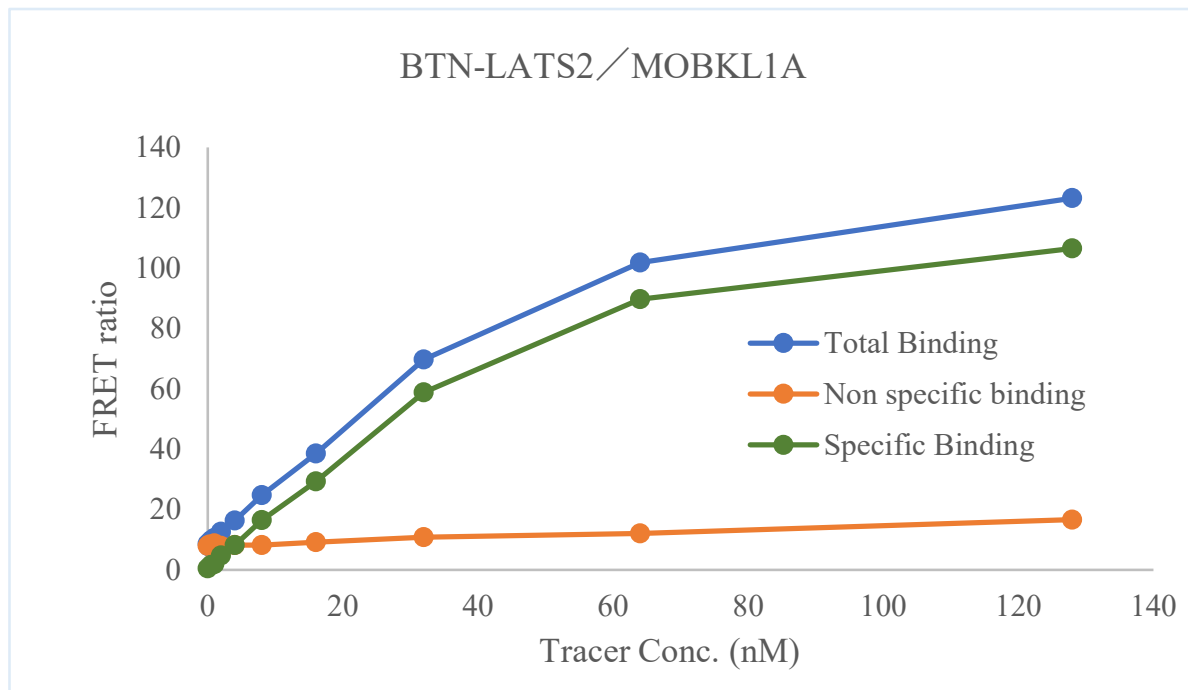


Kinase **BTN-LATS2/MOBKL1A**

Cat.# 01-424-20N

Tracer Revvity Inc.

Staurosporine-Red



KD (nM) 57.6

Kinase conc. 5nM

Eu labeled-streptavidin 5nM

Follow the guide to set up your HTRF® binding assay



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

