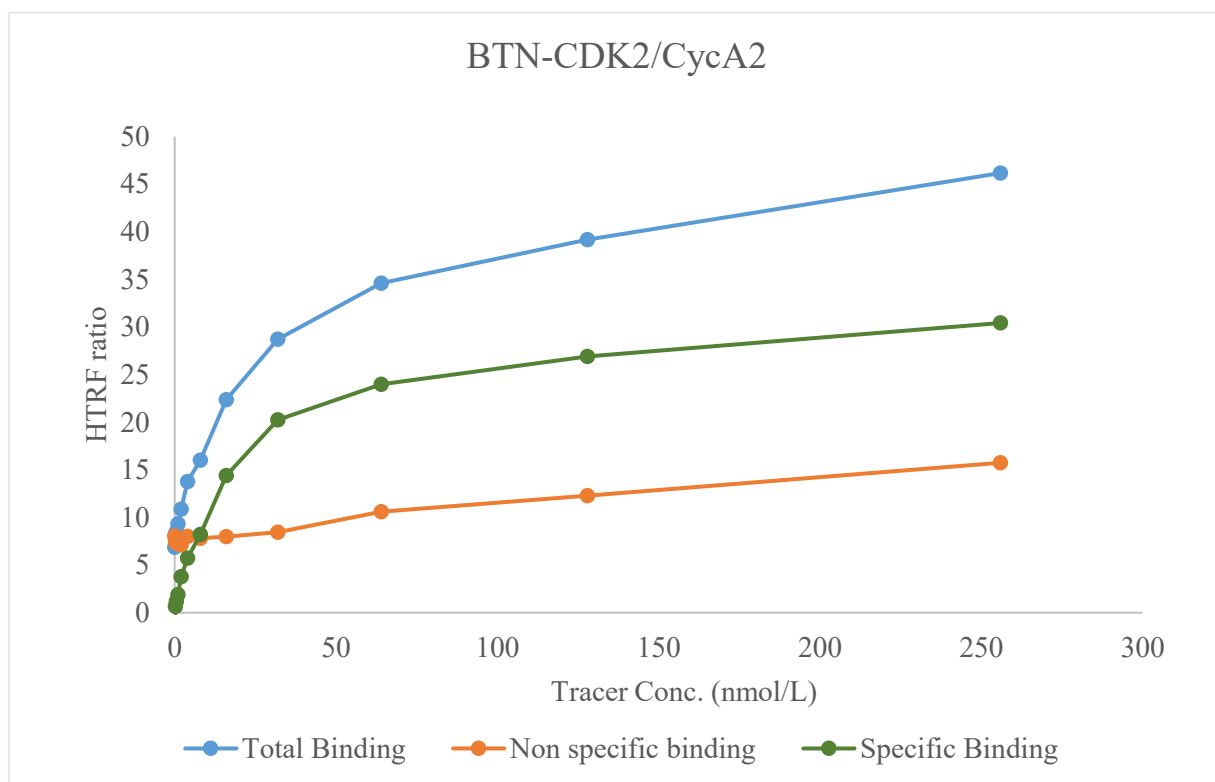


Kinase **BTN-CDK2/CycA2**

Cat.# 04-403-20N



Tracer

Thermo Scientific

Kinase Tracer 222, PV6121

KD

20 nmol/L

Kinase Conc.

5 nmol/L

Eu labeled-streptavidin

5 nmol/L

Metal

Mg

Metal Conc.

5 mmol/L

Additives

-

Reaction Time

1 Hour