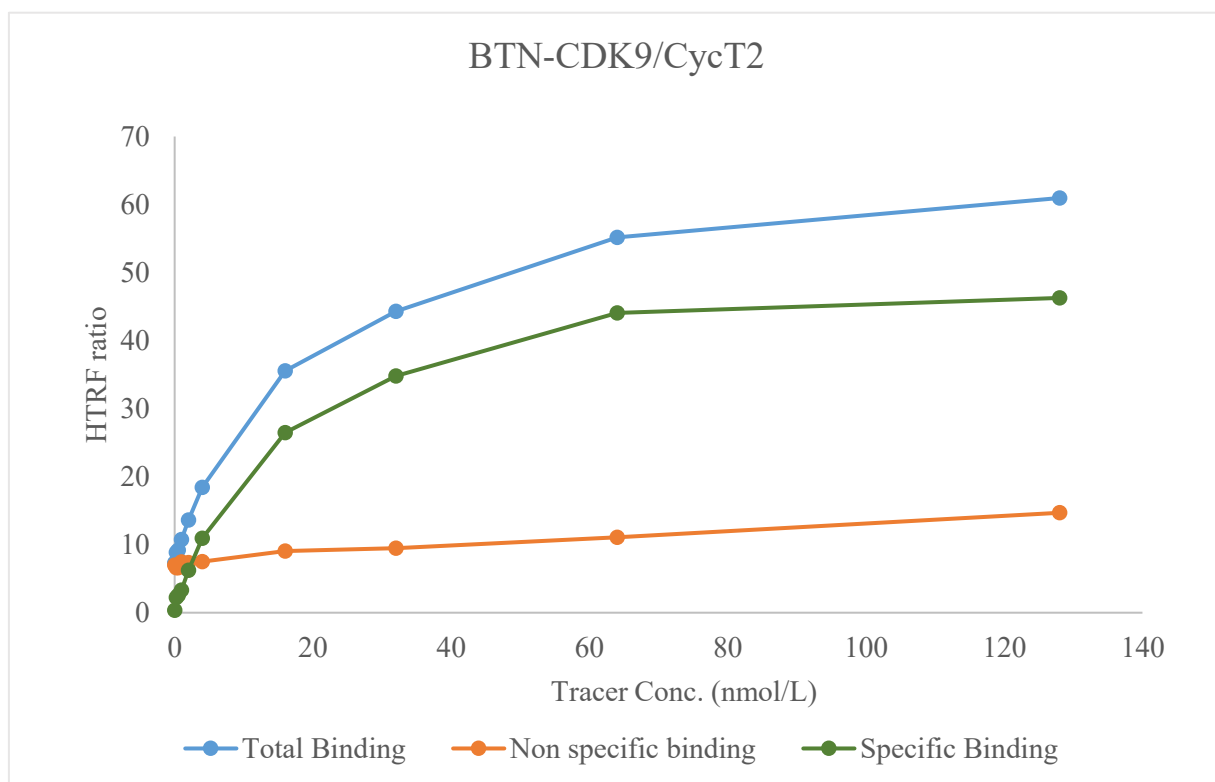


Kinase **BTN-CDK9/CycT2**

Cat.# 04-421-20N



Tracer

Thermo Scientific

Kinase Tracer 236, PV5592

KD

17 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