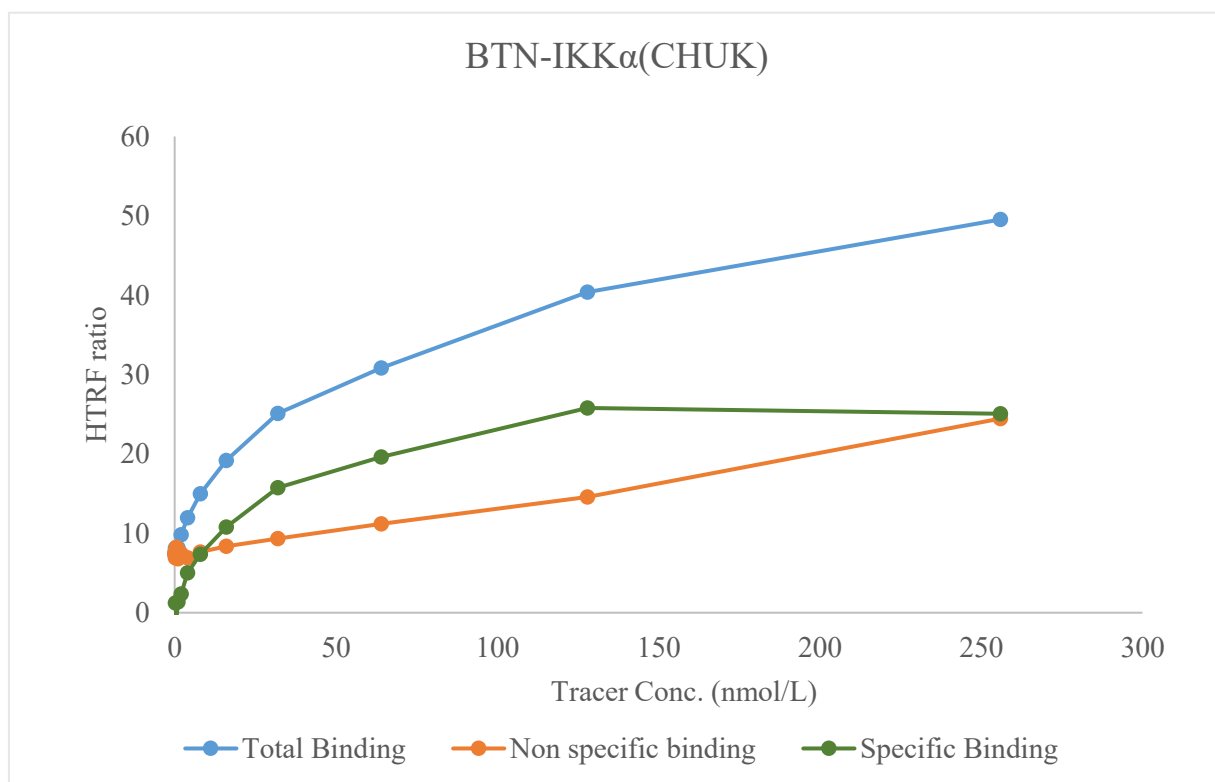


Kinase **BTN-IKK α (CHUK)**

Cat.# 05-412-20N



Tracer	Thermo Scientific
	Kinase Tracer 236, PV5592
KD	25 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