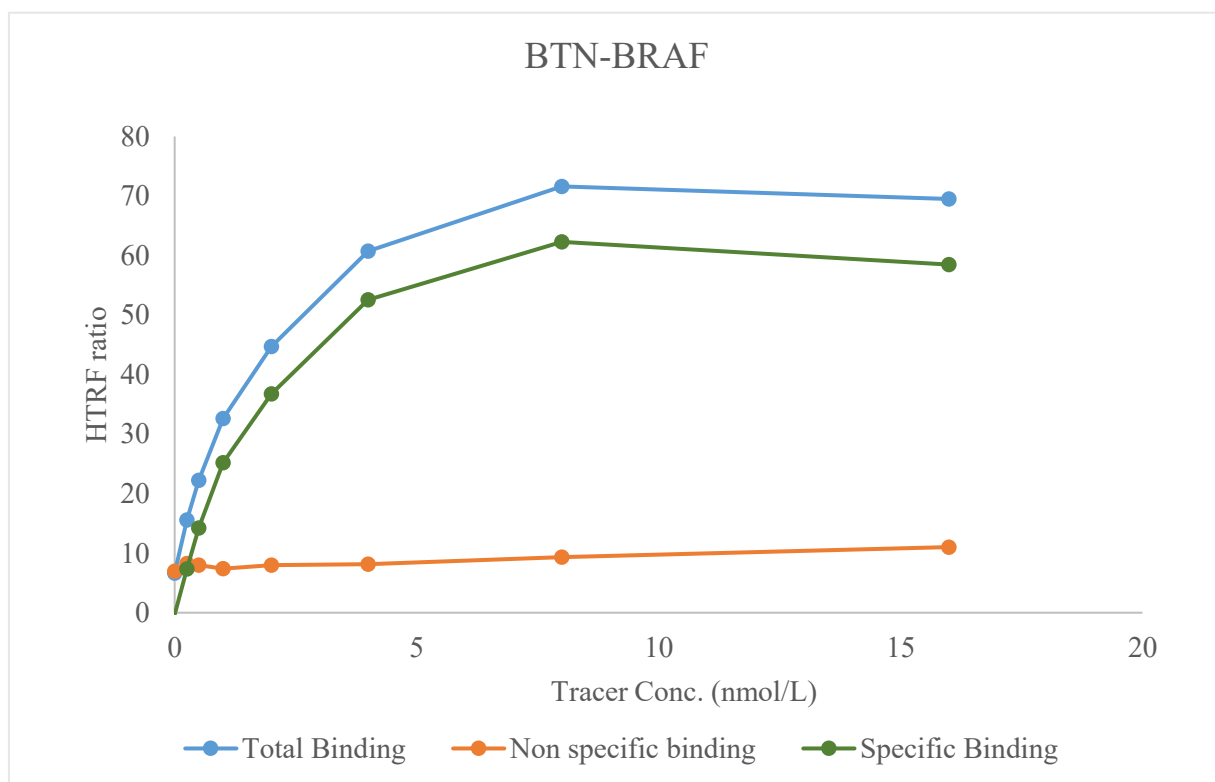


Kinase **BTN-BRAF**

Cat.# 09-422-20N



|                         |                           |
|-------------------------|---------------------------|
| Tracer                  | Thermo Scientific         |
|                         | Kinase Tracer 178, PV5593 |
| KD                      | 1.3 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                    |