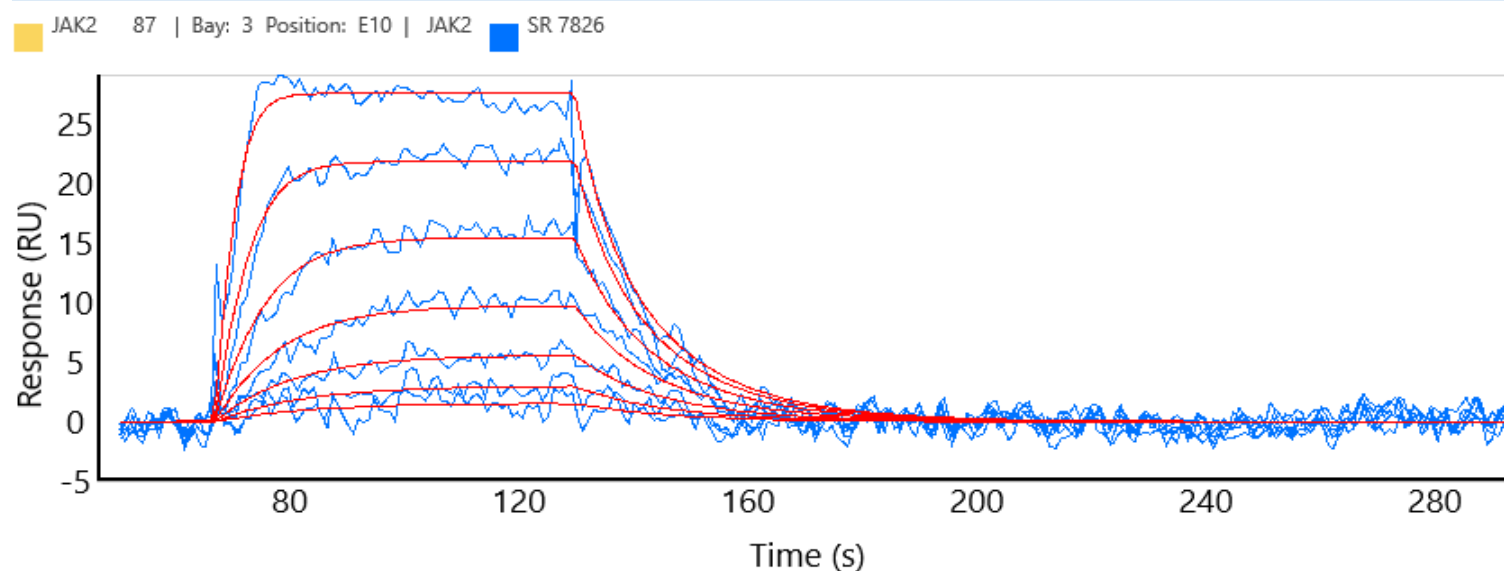


JAK2 binding kinetics on Carterra's HT-SPR platform

Product BTN-JAK2
Cat# 08-445-20N
Inhibitor SR 7826



Carterra's High Throughput SPR Platforms

The Carterra LSA^{XT} and Ultra utilize Surface Plasmon Resonance to detect binding interactions in real-time for up to 384 samples in parallel.

You can find more information at <https://carterra-bio.com/>

