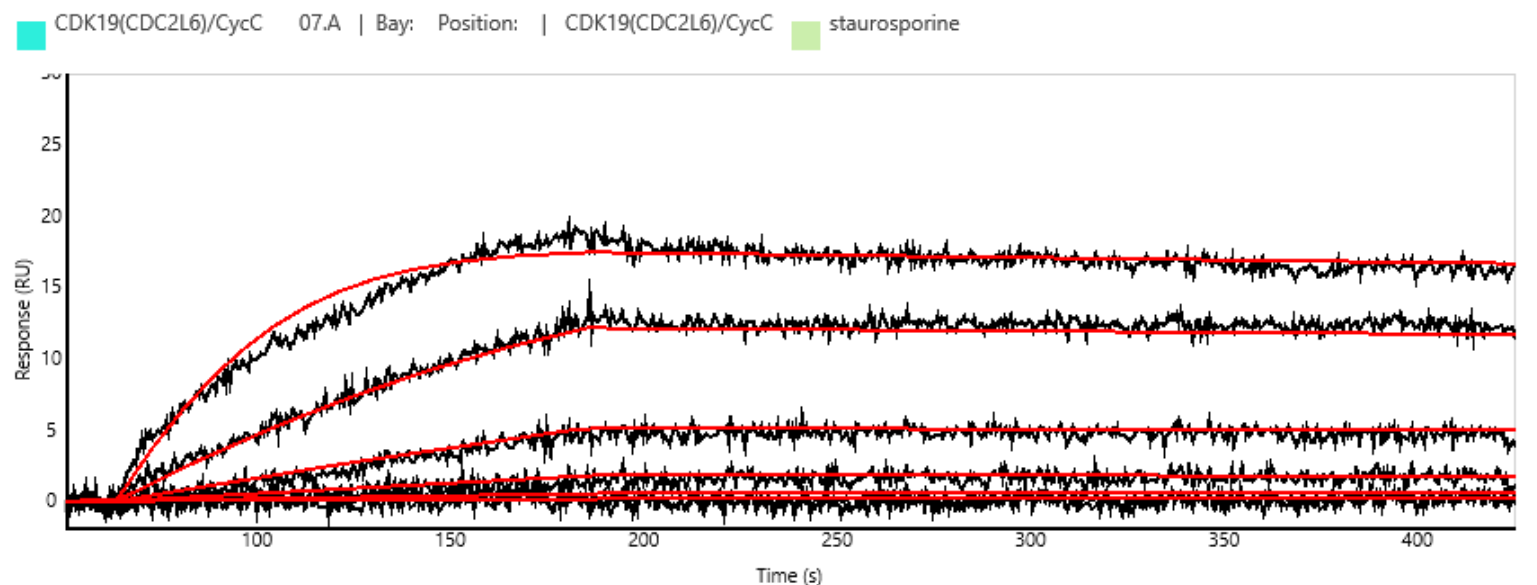


CDK19(CDC2L6)/CycC binding kinetics on Carterra's HT-SPR platform

Product BTN-CDK19(CDC2L6)/CycC

Cat# 04-412-20N

Inhibitor Staurosporine



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/>

