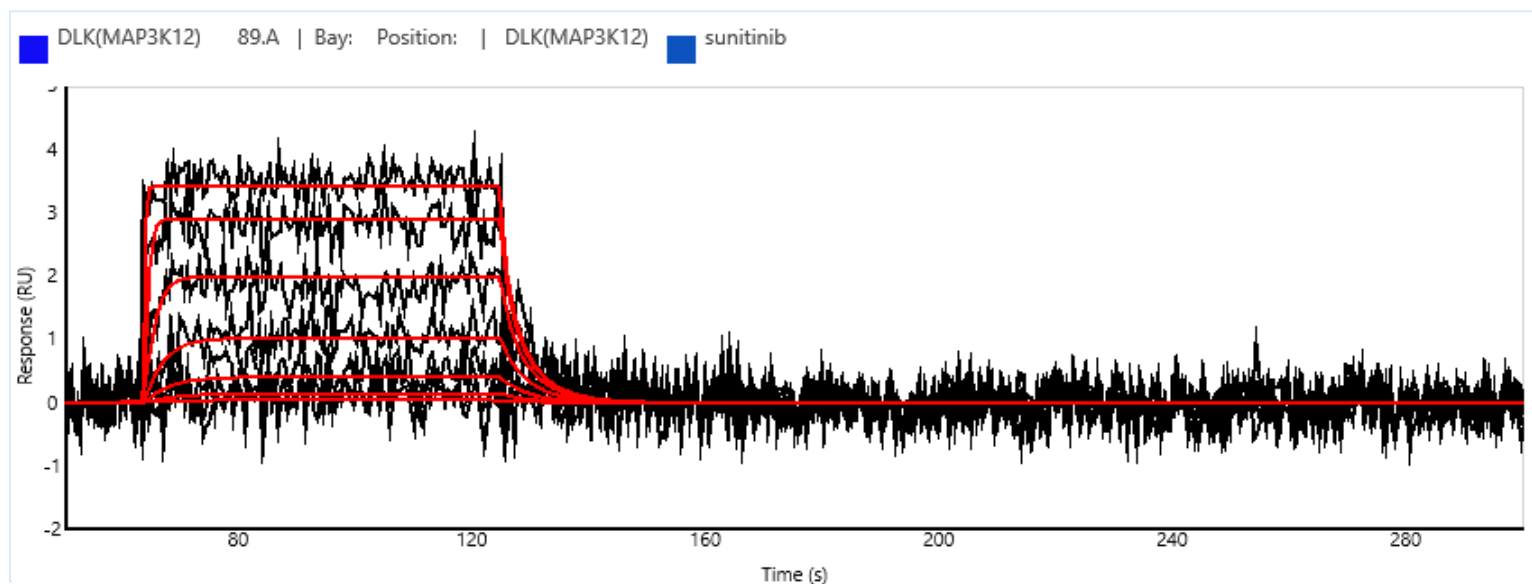


DLK(MAP3K12) binding kinetics on Carterra's HT-SPR platform

Product BTN-DLK(MAP3K12)

Cat# 09-411-20N

Inhibitor Sunitinib



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/>

