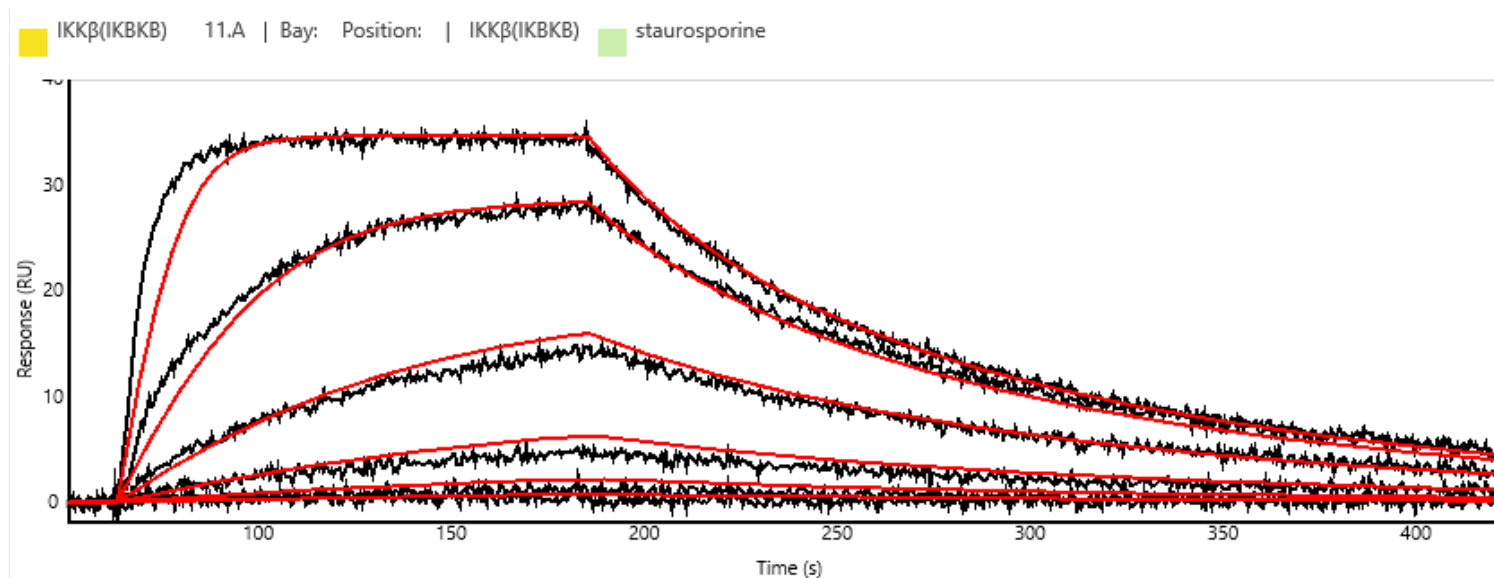


## IKK $\beta$ (IKBKB) binding kinetics on Carterra's HT-SPR platform

Product BTN-IKK  $\beta$  (IKBKB)

Cat# 05-484-20N

Inhibitor Staurosporine



## Carterra's High Throughput SPR Platforms

The Carterra LSA<sup>XT</sup> 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/>

