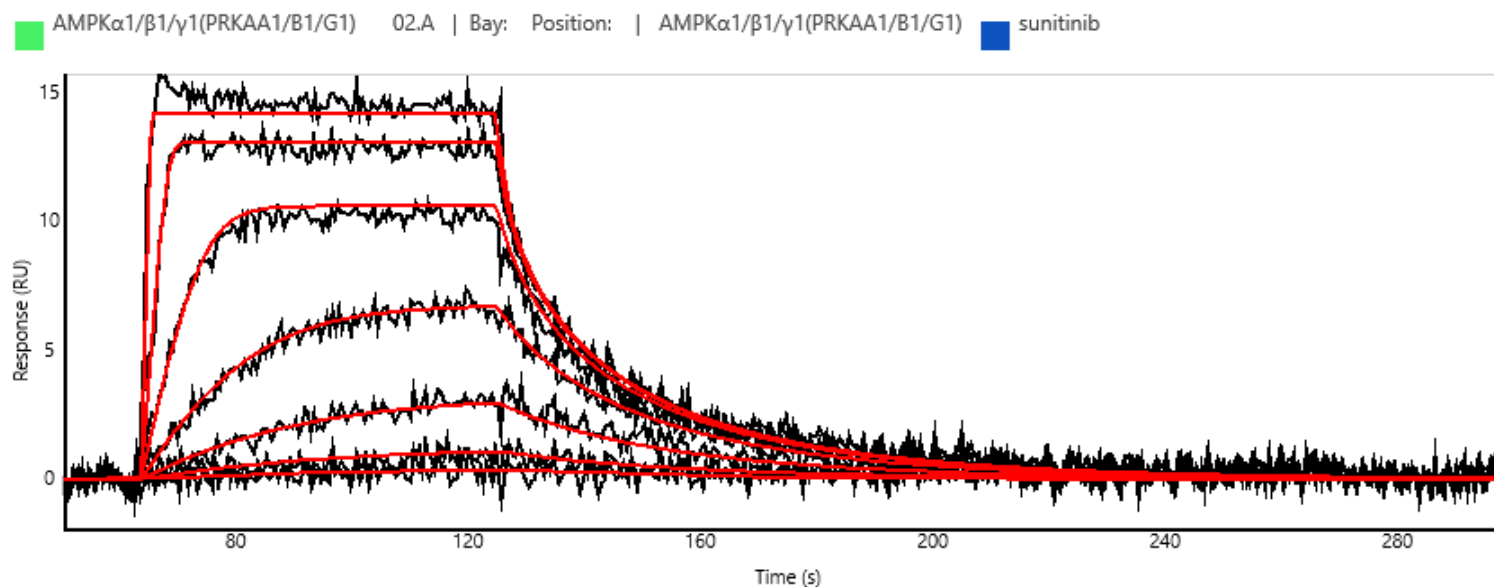


## AMPK $\alpha$ 1/ $\beta$ 1/ $\gamma$ 1 (PRKAA1/B1/G1) binding kinetics on Carterra's HT-SPR platform

Product BTN-AMPK $\alpha$ 1/ $\beta$ 1/ $\gamma$ 1 (PRKAA1/B1/G1)

Cat# 02-413-20N

Inhibitor Sunitinib



### 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/>

