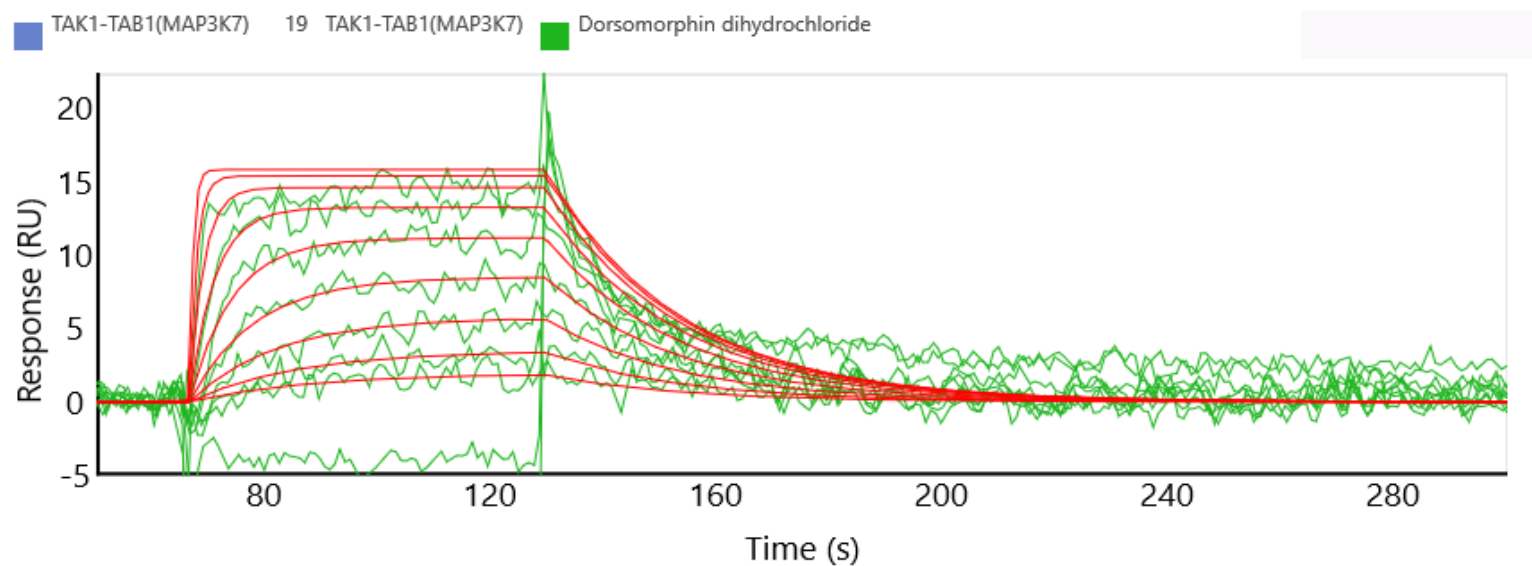


TAK1-TAB1 (MAP3K7) binding kinetics on Carterra's HT-SPR platform

Product BTN-TAK1-TAB1(MAP3K7)
 Cat# 09-419-21C
 Inhibitor Dorsomorphin dihydrochloride



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/>

