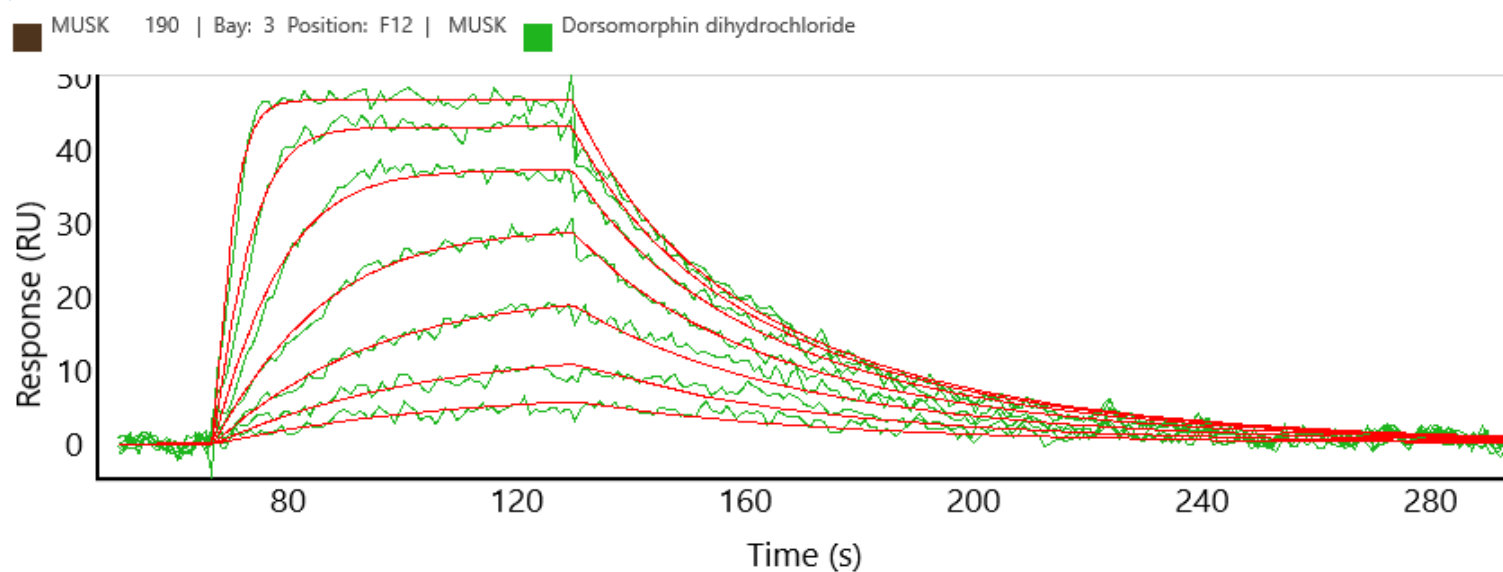


## MUSK binding kinetics on Carterra's HT-SPR platform

Product      BTN-MUSK  
Cat#          08-453-20N  
Inhibitor     Dorsomorphin dihydrochloride



## 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/>

