

**Sample name:** ADB-BUTINACA  
**Data file:** 29350-0576506.dx  
**Acq. method:** GC-MSD General Method 2.amx  
**Acquired on:** 2020-01-09 19:36:49-05:00  
**Method Description:**

**Operator :** Samantha Goodwin  
**Instrument:** ISO GCMS2  
**Inj. volume:** 1.000  
**Column:**

