

File :C:\ISO\23OCT2018\11151-0539796.D
Operator : JianmeiL
Acquired : 23 Oct 2018 16:10 using AcqMethod CAY_DRUG.M
Instrument : ISO_GCMS1
Sample Name: Mitragynine
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@240C, 30C/min to 300C
Vial Number: 8

