

File :C:\ISO\27JUL2017\22801-0511357.D
Operator : JianmeiL
Acquired : 27 Jul 2017 13:01 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: Crotonyl Fentanyl
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C
Vial Number: 7

