

File :C:\ISO\04OCT2018\17049-0541382.D
Operator : JianmeiL
Acquired : 4 Oct 2018 19:29 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: para-Fluorobutryl fentanyl (hydrochloride)
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C
Vial Number: 13

