

File :C:\ISO\Data\10JUN2015\14719-0467883.D
Operator : LauraN
Acquired : 10 Jun 2015 10:11 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: Fentanyl (HCl)
Misc Info : 30mx0.32mm,0.5uRtx-5MS,1min@50C-30C/min to 300C
Vial Number: 4

