

File :C:\ISO\12MAY2017\20207-0505593.D
Operator : JianmeiL
Acquired : 12 May 2017 18:45 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: ortho Fluoroisobutyl fentanyl (HCl)
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C
Vial Number: 9

