

File :C:\ISO\Data\24MAY2016\18705-0482873.D  
Operator : LauraN  
Acquired : 24 May 2016 14:06 using AcqMethod CAY\_DRUG\_CATH.M  
Instrument : ISO\_GCMS1  
Sample Name: Furanyl fentanyl (HCl)  
Misc Info : 30m x 0.32mm, Rtx-5MS, 1min@50C,30C/min to 300C  
Vial Number: 3

