

File :C:\ISO\16AUG2016\17049-0487736.D
Operator : NathanL
Acquired : 16 Aug 2016 13:58 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: para-Fluorobutyryl fentanyl (HCl)
Misc Info : 30m x 0.32mm, Rtx-5MS, 1min@50C,30C/min to 300C 1mg/ml in MeOH
Vial Number: 3

