

File :C:\ISO\15FEB2017\18089-0495945.D
Operator : LauraN
Acquired : 15 Feb 2017 18:16 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: 4-methoxy-Butyryl fentanyl (hydrochloride)
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C
Vial Number: 5

