

File :C:\ISO\21NOV2017-non ISO\23548-0516520R.D
Operator : JianmeiL
Acquired : 22 Nov 2017 12:13 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: ortho-methyl Acetyl fentanyl (hydrochloride)
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C
Vial Number: 10

