

File :C:\ISO\08AUG2017\22979-0512561.D
Operator : JianmeiL
Acquired : 8 Aug 2017 15:53 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: para-methyl Methoxyacetyl fentanyl (hydrochloride)
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C
Vial Number: 8

