

File :C:\ISO\29JAN2018-non ISO\14897-0521204 R.D  
Operator : JianmeiL  
Acquired : 31 Jan 2018 8:55 using AcqMethod CAY\_DRUG\_CATH.M  
Instrument : ISO\_GCMS1  
Sample Name: N-methyl-2-AI (hydrochloride)  
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C  
Vial Number: 18

