

File :C:\ISO\19APR2018-NON ISO\10643-0521253\_DRUG.D  
Operator : JianmeiL  
Acquired : 20 Apr 2018 12:23 using AcqMethod CAY\_DRUG.M  
Instrument : ISO\_GCMS1  
Sample Name: JWH 182  
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@240C,30C/min to 300C  
Vial Number: 16

