

File :C:\ISO\12OCT2017\15389-0507619_DRUG.D
Operator : JianmeiL
Acquired : 13 Oct 2017 8:44 using AcqMethod CAY_DRUG.M
Instrument : ISO_GCMS1
Sample Name: SDB-005
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@240C,30C/min to 300C
Vial Number: 3

