

File :C:\ISO\04OCT2018\90080-0542338.D
Operator : JianmeiL
Acquired : 4 Oct 2018 15:53 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: Cannabidiol
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C
Vial Number: 4

