

File :C:\ISO\26JAN2018\11488-0522537.D
Operator : JianmeiL
Acquired : 26 Jan 2018 15:55 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: Salvinorin B
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C
Vial Number: 4

