

File :C:\ISO\09OCT2018 ISO\16519-0541512.D
Operator : JianmeiL
Acquired : 9 Oct 2018 16:22 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: (R)-Ketamine HCl
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C
Vial Number: 4

