

File :C:\ISO\13JUN2018\10006350-0533475.D
Operator : JianmeiL
Acquired : 13 Jun 2018 22:51 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: HU-211
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C
Vial Number: 13

