

File :C:\ISO\26JUN2018\11565-0534002.D
Operator : JianmeiL
Acquired : 26 Jun 2018 13:01 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: XLR11
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C
Vial Number: 3

