

File :C:\ISO\29AUG2018 skb\16435-0530804.D
Operator : SBouhall
Acquired : 29 Aug 2018 11:33 using AcqMethod CAY_DRUG.M
Instrument : ISO_GCMS1
Sample Name: Flubromazolam
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@240C,30C/min to 300C
Vial Number: 3

