

File :C:\ISO\23JUL2018\20194-0533353.D
Operator : JianmeiL
Acquired : 23 Jul 2018 19:20 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: 4-cyano CUMYL-BUTINACA
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C
Vial Number: 15

