

File :C:\ISO\18OCT2016\14292-0490911.D
Operator : LauraN
Acquired : 19 Oct 2016 13:00 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: ADB-FUBINACA
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C
Vial Number: 14

