

File :C:\ISO\30MAR2017\14968-0496714_drug.D
Operator : SamanthaB
Acquired : 31 Mar 2017 13:28 using AcqMethod CAY_DRUG.M
Instrument : ISO_GCMS1
Sample Name: FDU-PB-22
Misc Info :
Vial Number: 3

