

File :Y:\Forensics\11312-0439953.D
Operator : PDK
Acquired : 30 Aug 2012 13:09 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: 4-hydroxy DiPT
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 17

