

File :Y:\Forensics\13885-0442837.D
Operator : PDK
Acquired : 4 Nov 2012 18:14 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: DOI
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 10

