

File :Y:\Forensics\14078-0451552.D
Operator : LJJ
Acquired : 13 Nov 2013 19:08 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: R-MMC, 3,4-Methylenedioxy-5-methyl-N-ethyl Ca
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 8

