

File :Y:\Forensics\13970-0444801-2.D
Operator : PDK
Acquired : 21 Mar 2013 14:30 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: 2,3-MDMA
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 11

