

File :Y:\Forensics\10007691-0438238-1.D
Operator : PDK
Acquired : 6 Jun 2012 15:26 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: URB754
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 11

