

File :Y:\Forensics\11048-0433283-2.D
Operator : PDK
Acquired : 3 Apr 2012 9:16 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: 2-AI
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 1

