

File :Y:\Forensics\11473-0436047.D
Operator : PDK
Acquired : 17 Aug 2012 12:25 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: 2C-B-fly
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 9

