

File :Y:\Forensics\11008-0436043.D
Operator : PDK
Acquired : 4 Apr 2012 10:40 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: AM2233
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 3

