

File :Y:\Forensics\10567-0439464.D
Operator : PDK
Acquired : 8 Aug 2012 19:55 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: AM694
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 12

