

File :Y:\Forensics\10707-0430471.D
Operator : PDK
Acquired : 20 Aug 2012 20:17 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: AM2201
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 17

