

File :Y:\Forensics\11282-0448969.D
Operator : PDK
Acquired : 16 Aug 2013 23:02 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: AM1248
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 6

