

File :Y:\Forensics\11564-0446346.D
Operator : PDK
Acquired : 30 Apr 2013 21:05 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: STS-135
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 1

