

File :Y:\Forensics\11556-0441294.D
Operator : PDK
Acquired : 7 Nov 2012 13:15 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: N,N-Dimethylamphetamine
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 7

