

File :Y:\Forensics\11562-0438332-2.D
Operator : PDK
Acquired : 3 Aug 2012 20:44 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: para-Methoxymethamphetamine
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 17

