

File :Y:\Forensics\11139-0440299.D
Operator : PDK
Acquired : 17 Aug 2012 12:00 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: Methoxetamine
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-300C/min-300C
Vial Number: 8

