

File :Y:\Forensics\11200-0440095.D
Operator : PDK
Acquired : 22 Aug 2012 13:30 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: 2C-T-7
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-300C/min-300C
Vial Number: 2

