

File :Y:\Forensics\9001070-0439381-4.D
Operator : PDK
Acquired : 3 Aug 2012 16:33 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: 4-Fluoromethamphetamine
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 6

