

File :Y:\Forensics\9001863-0453399-2.D
Operator : PDK
Acquired : 11 Apr 2014 14:04 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: 2-Ethylamino-1-phenylbutane
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 5

