

File :Y:\Forensics\9001037-0440593.D
Operator : PDK
Acquired : 4 Nov 2012 23:54 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: JWH 122 8-methylnaphthyl isomer
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 3

