

File :Y:\Forensics\10828-0441500.D
Operator : PDK
Acquired : 5 Nov 2012 2:11 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: JWH 369
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 7

