

File :Y:\FORENSICS\13169-0437763.D
Operator : PDK
Acquired : 20 Apr 2012 19:04 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: JWH 018
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 9

