

File :Y:\Forensics\10009018-0437213.D
Operator : PDK
Acquired : 5 Nov 2012 00:29 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: JWH 015
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 5

