

File :Y:\Forensics\11201-0448879-2.D
Operator : PDK
Acquired : 17 Jul 2013 16:48 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: JWH 175
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 4

