

File :Y:\Forensics\9001002-0449146-3.D
Operator : PDK
Acquired : 20 Aug 2013 15:17 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: JWH 018 N-(1-methylbutyl) isomer
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 3

