

File :Y:\Forensics\RMS-1211-122-15-2.D
Operator : PDK
Acquired : 14 Oct 2013 16:32 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: JWH 018 benzimidazole analog
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 2

