

File :Y:\Forensics\10521-0443464.D
Operator : PDK
Acquired : 17 Dec 2012 22:53 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: JWH 018 N-(5-chloropentyl) analog
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 17

