

File :Y:\Forensics\10591-0432767.D
Operator : PDK
Acquired : 17 Sep 2012 19:42 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: JWH 122
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 11

