

File :Y:\Forensics\14901-0455692.D
Operator : PDK
Acquired : 15 Apr 2014 18:45 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: Pyrazolam
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 10

