

File :Y:\Forensics\15536-0456474.D
Operator : PDK
Acquired : 17 Apr 2014 00:13 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: 5-fluoro NPB-22
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 12

