

File :Y:\Forensics\14816-0452997.D
Operator : LJJ
Acquired : 14 Nov 2013 19:44 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: 5-fluoro MN-18
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 11

