

File :Y:\Forensics\15202-0453507.D  
Operator : LJJ  
Acquired : 22 Nov 2013 12:26 using AcqMethod CAY\_DRUG.M  
Instrument : Instrument #1  
Sample Name: AM2201 benzimidazole analog  
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C  
Vial Number: 18

