

File :Y:\Forensics\14789-0457719.D  
Operator : PDK  
Acquired : 24 Jun 2014 14:13 using AcqMethod CAY\_DRUG.M  
Instrument : Instrument #1  
Sample Name: THJ 2201  
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C  
Vial Number: 1

