

File :Y:\Forensics\14196-0450293-3.D  
Operator : PDK  
Acquired : 11 Sep 2013 15:41 using AcqMethod CAY\_DRUG\_CATH.M  
Instrument : Instrument #1  
Sample Name: HMA  
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C  
Vial Number: 16

