

File :Y:\Forensics\12041-0441193.D  
Operator : PDK  
Acquired : 19 Sep 2012 13:26 using AcqMethod CAY\_DRUG\_CATH.M  
Instrument : Instrument #1  
Sample Name: para-Methoxyamphetamine  
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C  
Vial Number: 13

