

File :Y:\Forensics\14084-0445105-3.D  
Operator : PDK  
Acquired : 25 Mar 2013 20:20 using AcqMethod CAY\_DRUG.M  
Instrument : Instrument #1  
Sample Name: JWH 073 6-methoxyindole analog  
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C  
Vial Number: 15

