

File :Y:\Forensics\11242-0443093-2.D
Operator : PDK
Acquired : 19 Dec 2012 16:24 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: 2,4-Dimethylethcathinone
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 9

