

File :Y:\Forensics\11243-0443056.D
Operator : PDK
Acquired : 19 Dec 2012 15:10 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: 2,4-Dimethylmethcathinone
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 8

