

File :Y:\Agilent GCMS\Forensics\11207-0434033.D  
Operator : PDK  
Acquired : 19 Mar 2012 15:04 using AcqMethod CAY\_DRUG\_CATH.M  
Instrument : Instrument #1  
Sample Name: 4-ethyl-N,N-Dimethylcathinone  
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C  
Vial Number: 5

