

File :Y:\Agilent GCMS\Forensics\11221-0436091.D
Operator : PDK
Acquired : 15 Mar 2012 13:34 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: 2-Methylethcathinone
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 6

