

File :Y:\Agilent GCMS\Forensics\9001034-0431263.D
Operator : PDK
Acquired : 21 Mar 2012 22:09 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: JWH 122 5-methylnaphthyl isomer
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 12

