

File :Y:\Agilent GCMS\Forensics\9001013-0434161.D
Operator : PDK
Acquired : 20 Mar 2012 12:36 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: JWH-073 N-(1,1-dimethylethyl) isomer
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 5

