

File :Y:\Agilent GCMS\Forensics\9001044-0431076.D  
Operator : PDK  
Acquired : 21 Mar 2012 23:50 using AcqMethod CAY\_DRUG.M  
Instrument : Instrument #1  
Sample Name: JWH 081 2-methoxynaphthyl isomer  
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C  
Vial Number: 15

