

File :Y:\Agilent GCMS\Forensics\9001045-0431272.D
Operator : PDK
Acquired : 23 Mar 2012 16:49 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: JWH 081 3-methoxynaphthyl isomer
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 19

