

File :Y:\Agilent GCMS\Forensics\10691-0427677.D
Operator : PDK
Acquired : 29 Feb 2012 20:44 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: JWH 018-N-(3-methylbutyl) isomer
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 10

