

File :Y:\Agilent GCMS\Forensics\13633-0416401.D
Operator : PDK
Acquired : 14 Feb 2012 16:39 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: JWH-019
Misc Info : 30mx0.32mm,0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 5

