

File :Y:\Agilent GCMS\Forensics\9001076-0430975.D
Operator : PDK
Acquired : 26 Mar 2012 10:28 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: JWH 073 4-methoxynaphthyl analog
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 4

