

File :Y:\Agilent GCMS\Forensics\9001004-0429777.D
Operator : PDK
Acquired : 21 Feb 2012 10:23 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: JWH 018 2'-naphthyl isomer
Misc Info : 30mx0.32mm,0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 1

