

File :Y:\Agilent GCMS\Forensics\10865-0429305.D
Operator : PDK
Acquired : 16 Feb 2012 23:19 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: RCS-4 2-methoxy isomer
Misc Info : 30mx0.32mm,0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 11

