

File :Y:\Agilent GCMS\Forensics\10798-0427940.D
Operator : PDK
Acquired : 16 Mar 2012 12:44 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: RCS-4-C4-homolog
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 5

