

File :Y:\Agilent GCMS\Forensics\10645-0424539.D
Operator : PDK
Acquired : 1 Mar 2012 20:46 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: RCS-4
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 7

