

File :Y:\Agilent GCMS\Forensics\10869-0429225-3.D
Operator : PDK
Acquired : 19 Mar 2012 12:02 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: AM694 4-iodo isomer
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 7

