

File :Y:\Agilent GCMS\Forensics\11206-0435856.D
Operator : PDK
Acquired : 19 Mar 2012 20:59 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: 1,4-DibenZyl piperazine
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 14

