

File :Y:\Agilent GCMS\Forensics\10918-0428460-2.D
Operator : PDK
Acquired : 14 Mar 2012 13:23 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: (±)3-epi CP 47,497
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 9

