

File :Y:\Agilent GCMS\Forensics\10889-0430308.D
Operator : PDK
Acquired : 16 Mar 2012 18:13 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: CP 47,497-para-quinone analog
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 19

