

File :Y:\Agilent GCMS\Forensics\9001187-0435064.D
Operator : PDK
Acquired : 23 Mar 2012 19:35 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: 3-Methoxymethcathinone
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 11

