

File :Y:\Agilent GCMS\Forensics\10644-0425787.D
Operator : PDK
Acquired : 14 Feb 2012 20:02 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: JWH-210
Misc Info : 30mx0.32mm,0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 12

