

File :Y:\Agilent GCMS\Forensics\10697-0425868.D
Operator : PDK
Acquired : 12 Mar 2012 14:42 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: JWH 018 6-methoxyindole analog
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 7

