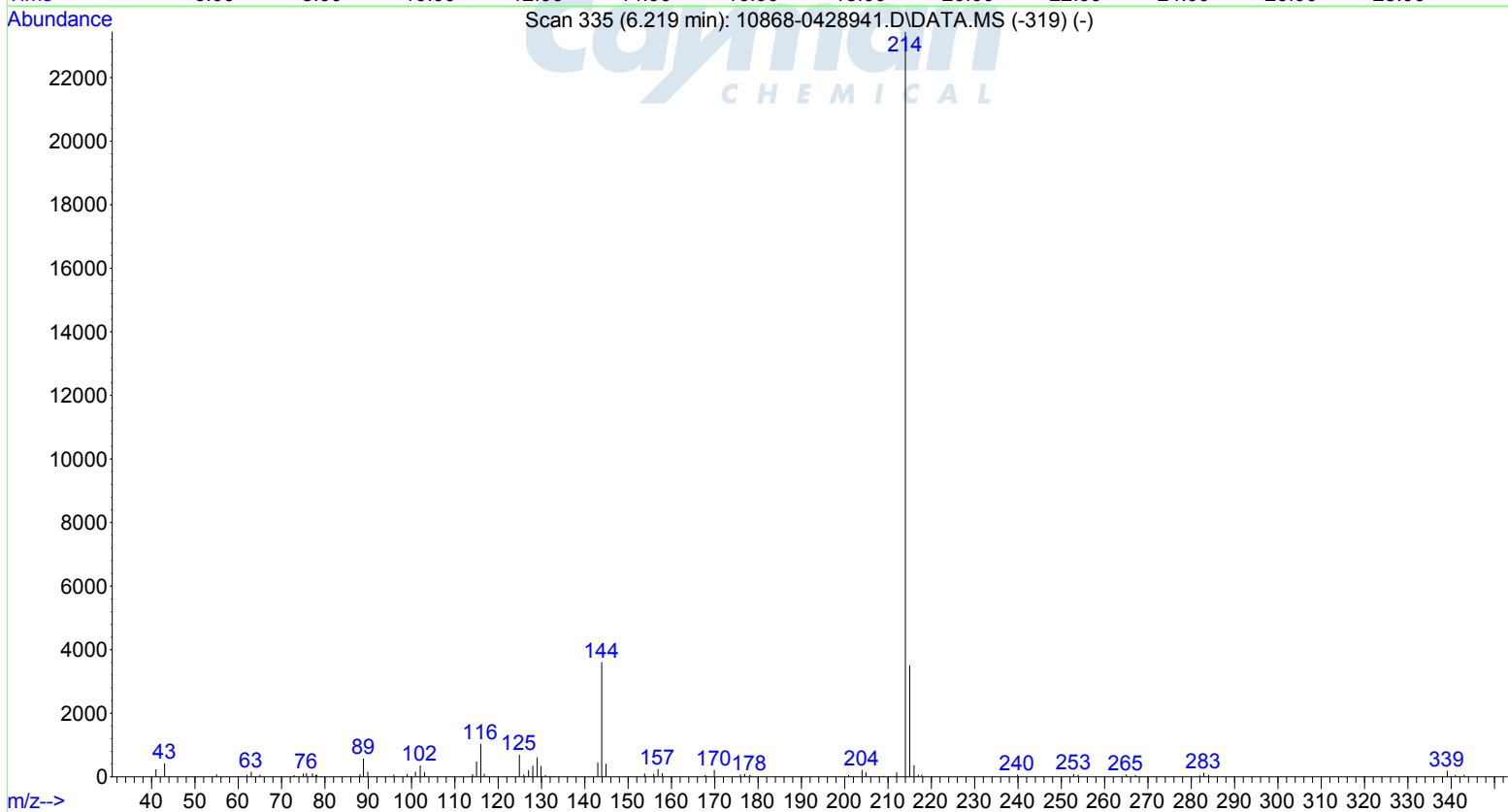
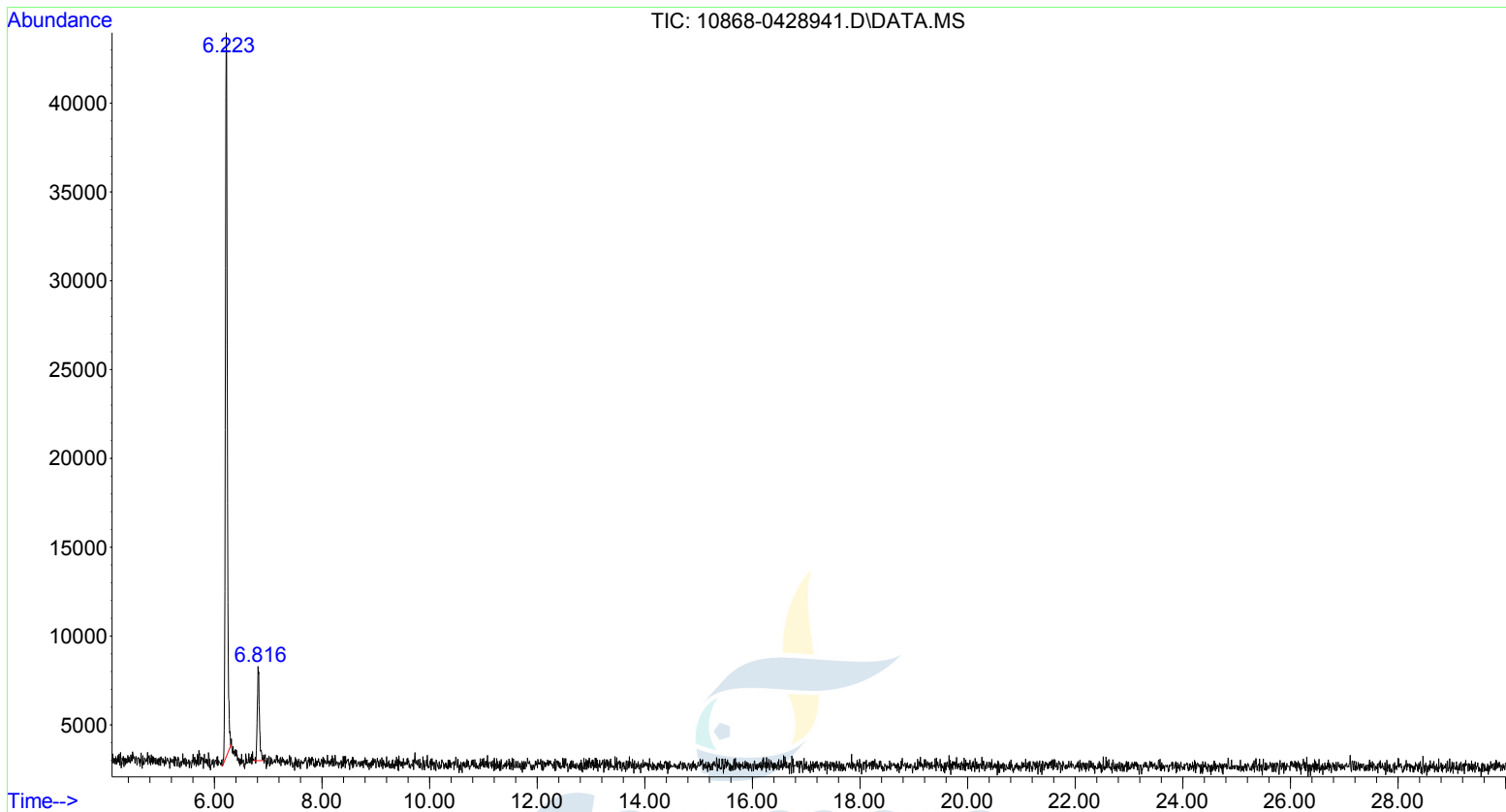


File :Y:\Agilent GCMS\Forensics\10868-0428941.D  
Operator : PDK  
Acquired : 16 Feb 2012 18:14 using AcqMethod CAY\_DRUG.M  
Instrument : Instrument #1  
Sample Name: JWH 203 3-chlorophenyl isomer  
Misc Info : 30mx0.32mm,0.5u Rtx-5MS, 240C-30C/min-300C  
Vial Number: 2



File :Y:\Agilent GCMS\Forensics\10868-0428941.D  
Operator : PDK  
Acquired : 16 Feb 2012 18:14 using AcqMethod CAY\_DRUG.M  
Instrument : Instrument #1  
Sample Name: JWH 203 3-chlorophenyl isomer  
Misc Info : 30mx0.32mm,0.5u Rtx-5MS, 240C-30C/min-300C  
Vial Number: 2

