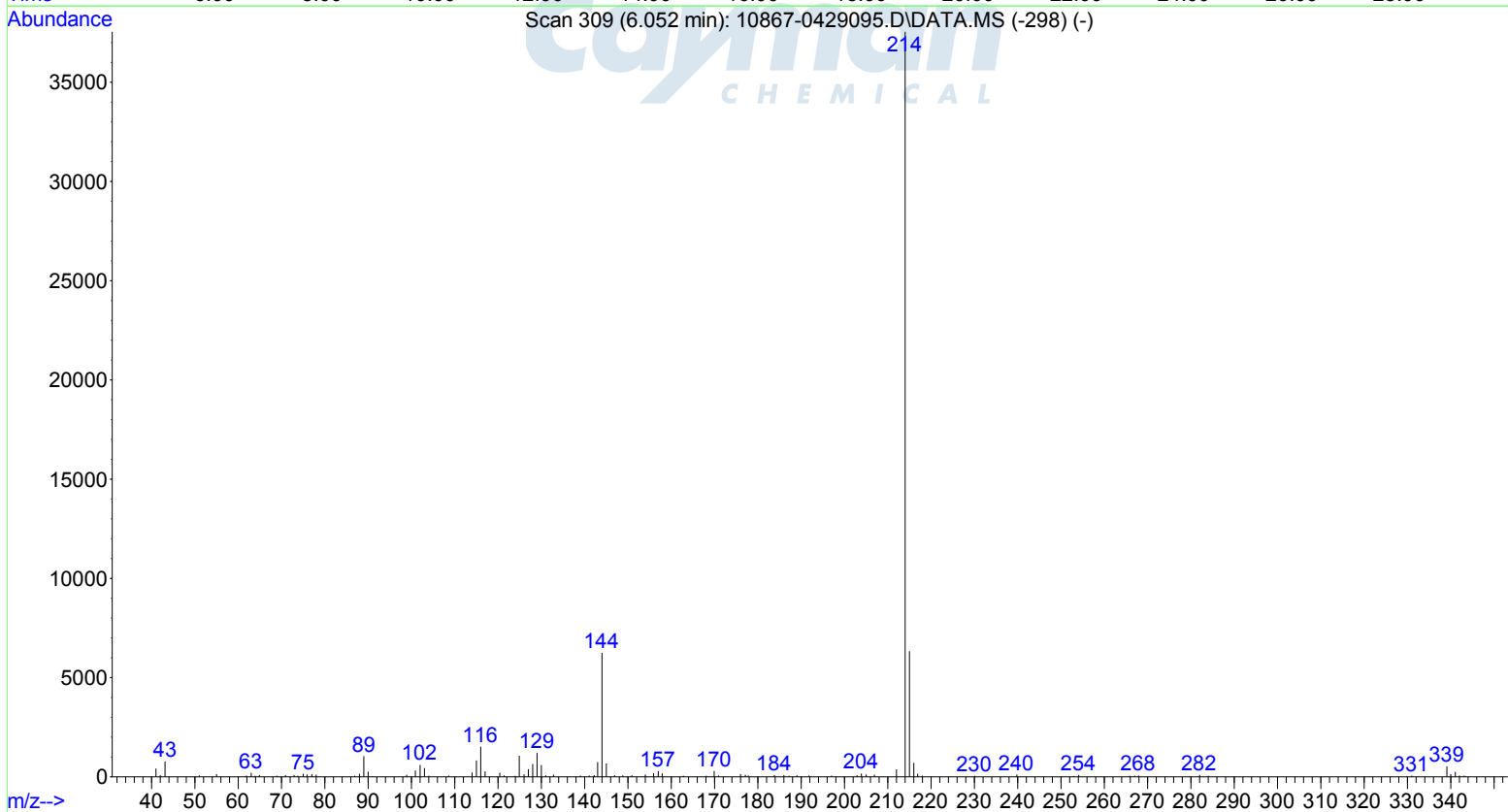
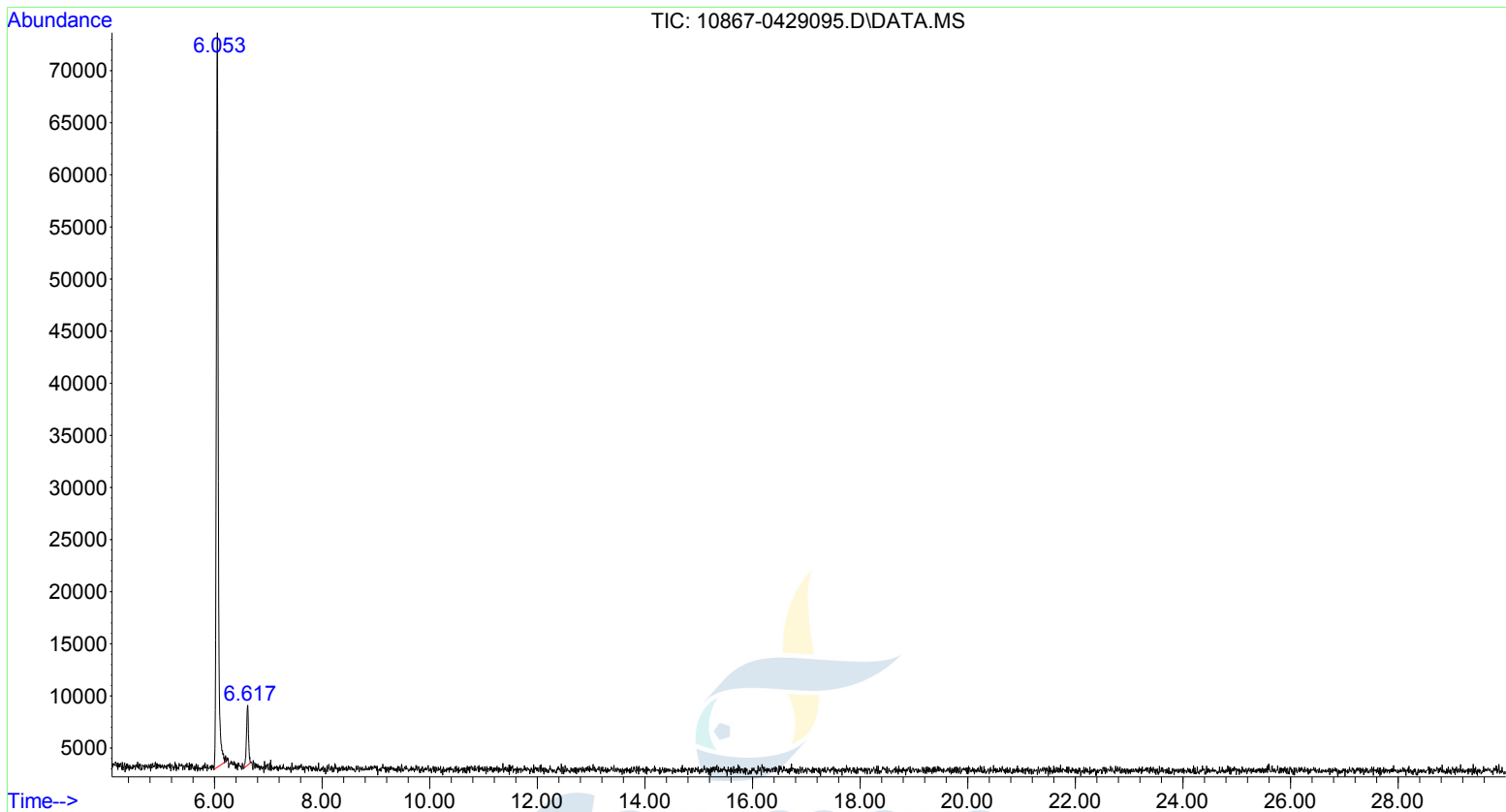


File :Y:\Agilent GCMS\Forensics\10867-0429095.D
Operator : PDK
Acquired : 16 Feb 2012 19:22 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: JWH 203 4-chlorophenyl isomer
Misc Info : 30mx0.32mm,0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 4



File :Y:\Agilent GCMS\Forensics\10867-0429095.D
Operator : PDK
Acquired : 16 Feb 2012 19:22 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: JWH 203 4-chlorophenyl isomer
Misc Info : 30mx0.32mm,0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 4

