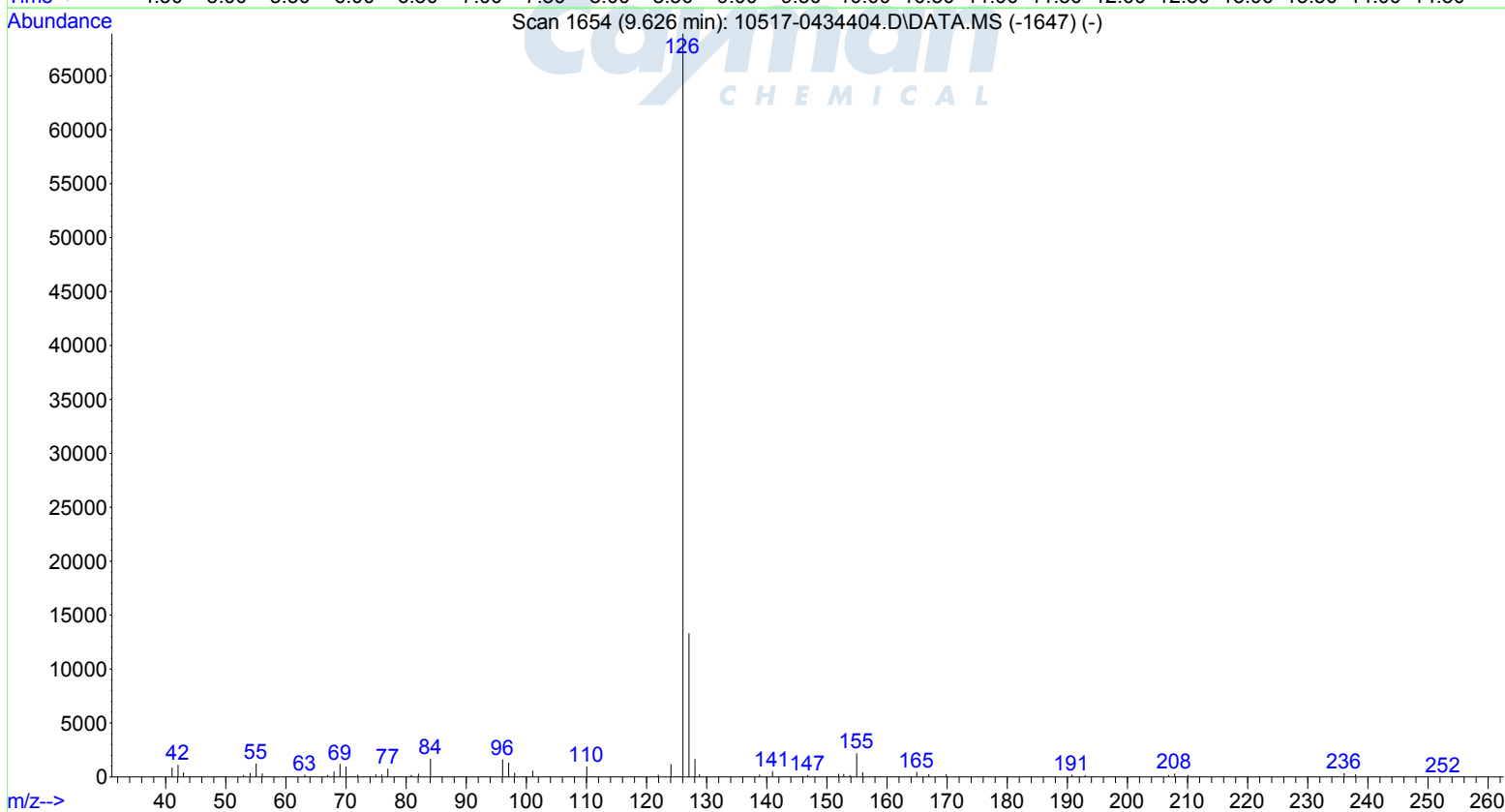
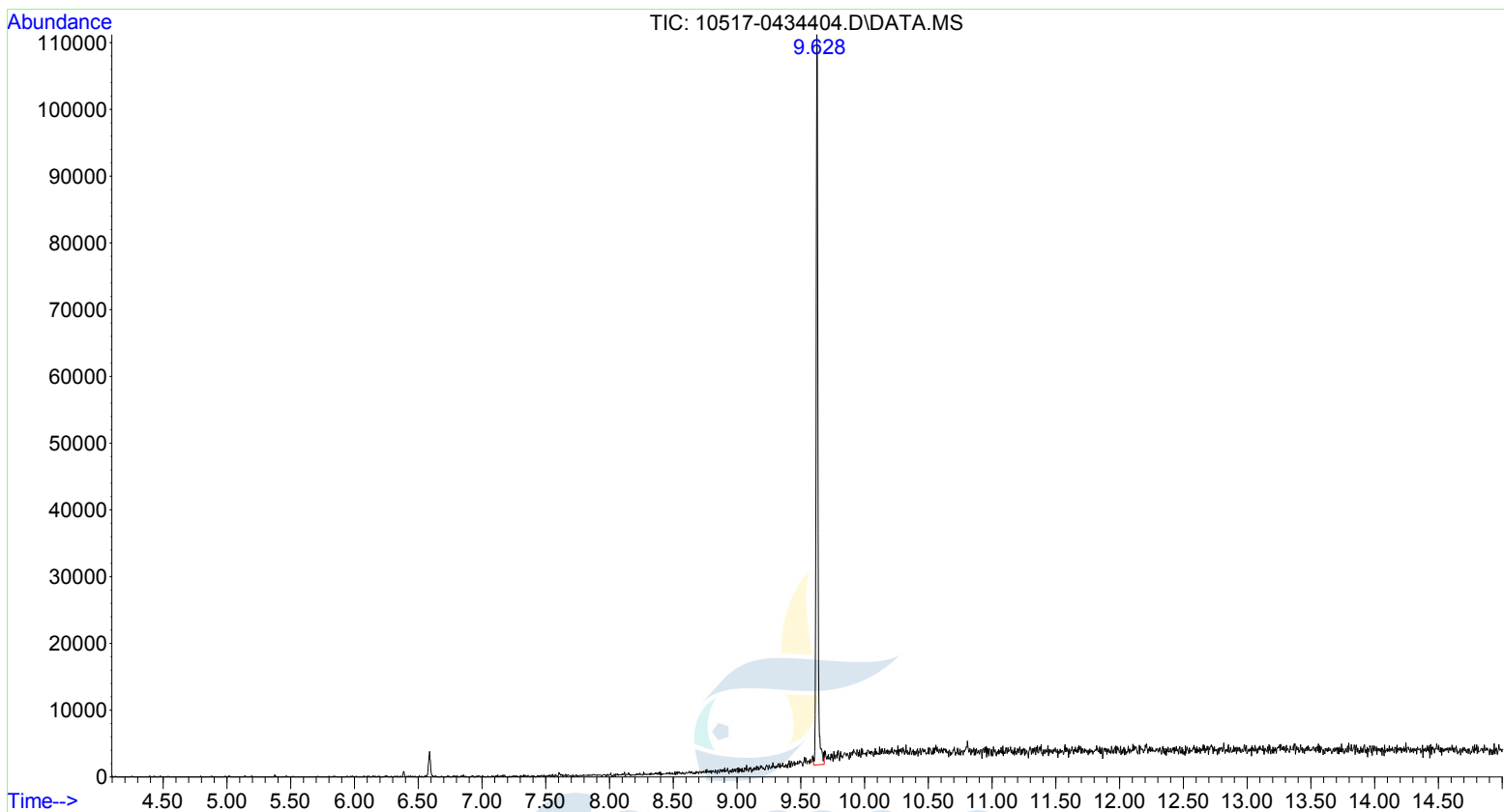


File :Y:\Agilent GCMS\Forensics\10517-0434404.D
Operator : PDK
Acquired : 29 Feb 2012 17:11 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: Naphyrone (HCl)
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 3



File :Y:\Agilent GCMS\Forensics\10517-0434404.D
Operator : PDK
Acquired : 29 Feb 2012 17:11 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: Naphyrone (HCl)
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 3

