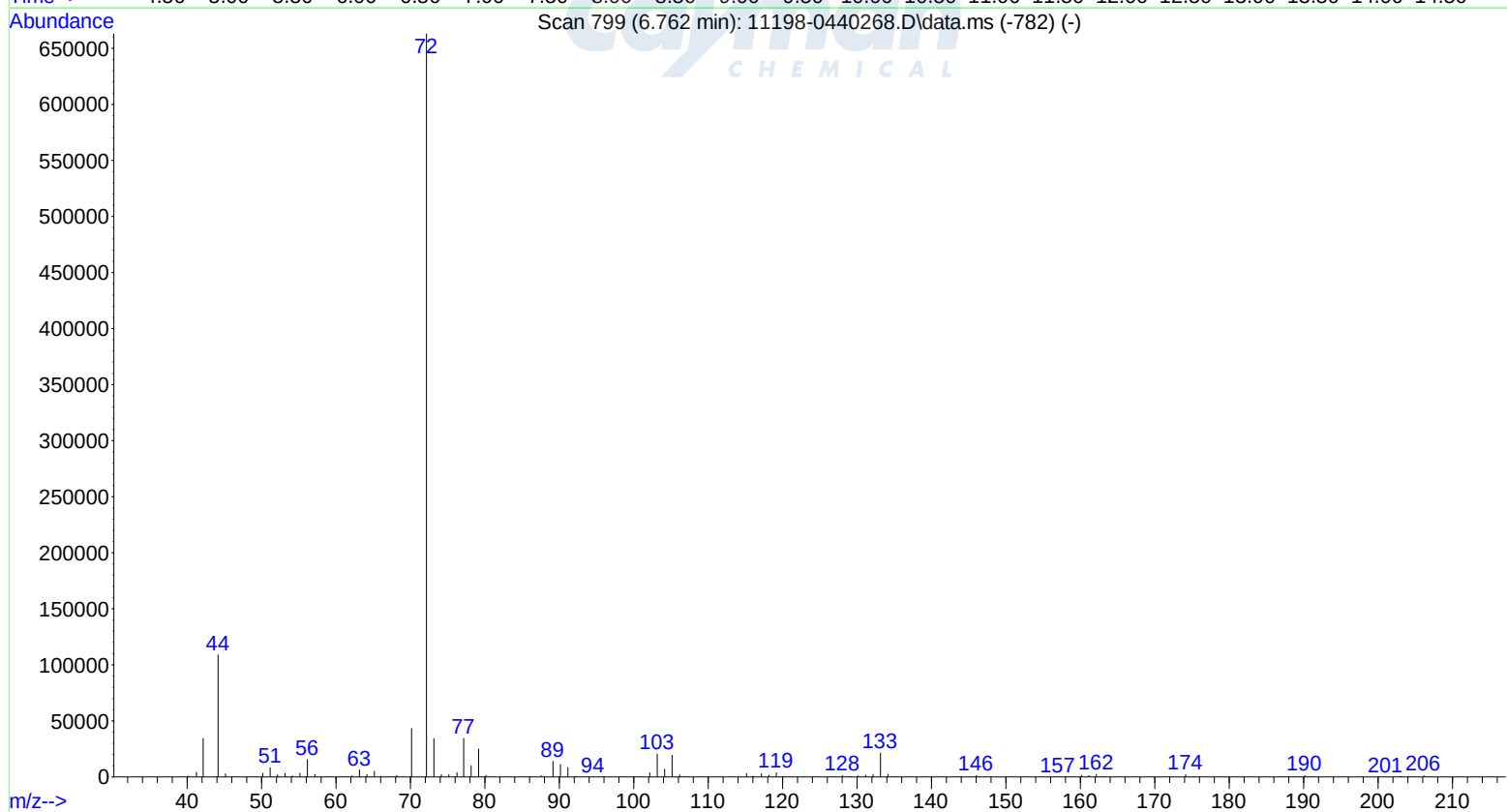
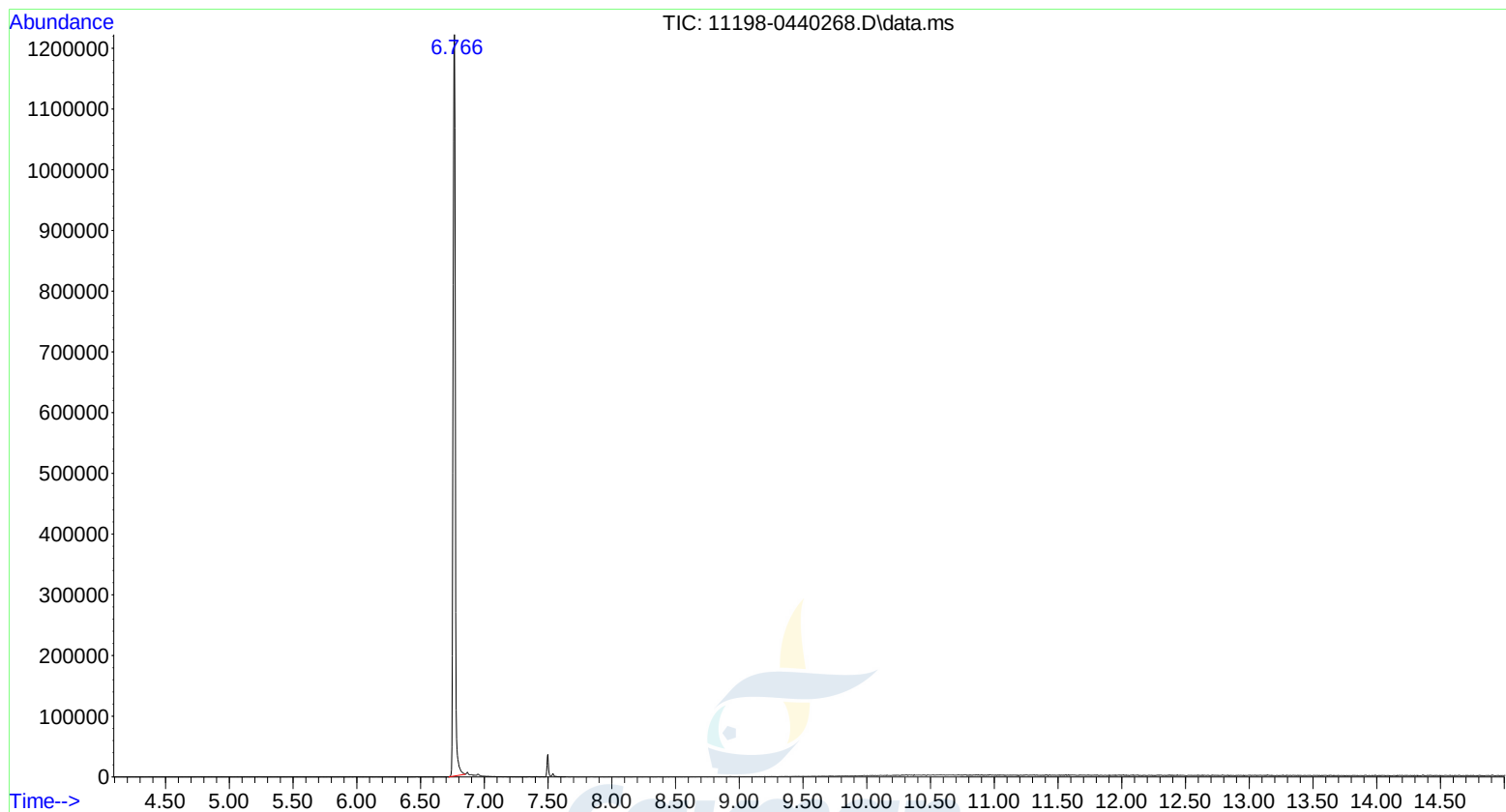


File :Y:\Forensics\11198-0440268.D
Operator : PDK
Acquired : 29 Aug 2012 15:58 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: 3-Ethylethcathinone
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 5



File :Y:\Forensics\11198-0440268.D

Operator : PDK

Acquired : 29 Aug 2012 15:58 using AcqMethod CAY_DRUG_CATH.M

Instrument : Instrument #1

Sample Name: 3-Ethylethcathinone

Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C

Vial Number: 5

