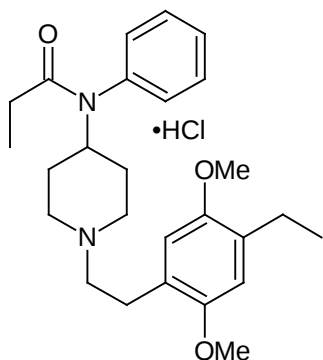


BATTELLE CERTIFICATE OF ANALYSIS



Name: N-(2C-E) fentanyl HCl

Supplier: Cayman Chemical

Catalog No.: 28973

Lot #: N/A

Chemical Formula: $C_{26}H_{36}N_2O_3$ (free base)
 $C_{26}H_{37}ClN_2O_3$ (salt)

Molecular Weight: 424.58 (free base)
461.04 (salt)

Analytical Results

Monoisotopic Exact 425.2799 (calc. for $[M+H]^+$)

Mass: 425.2798 (obsd. ESI-MS)

Isotopic Purity: N/A

Date of COA: April 21, 2020

COA Prepared by: E. A. Lucas

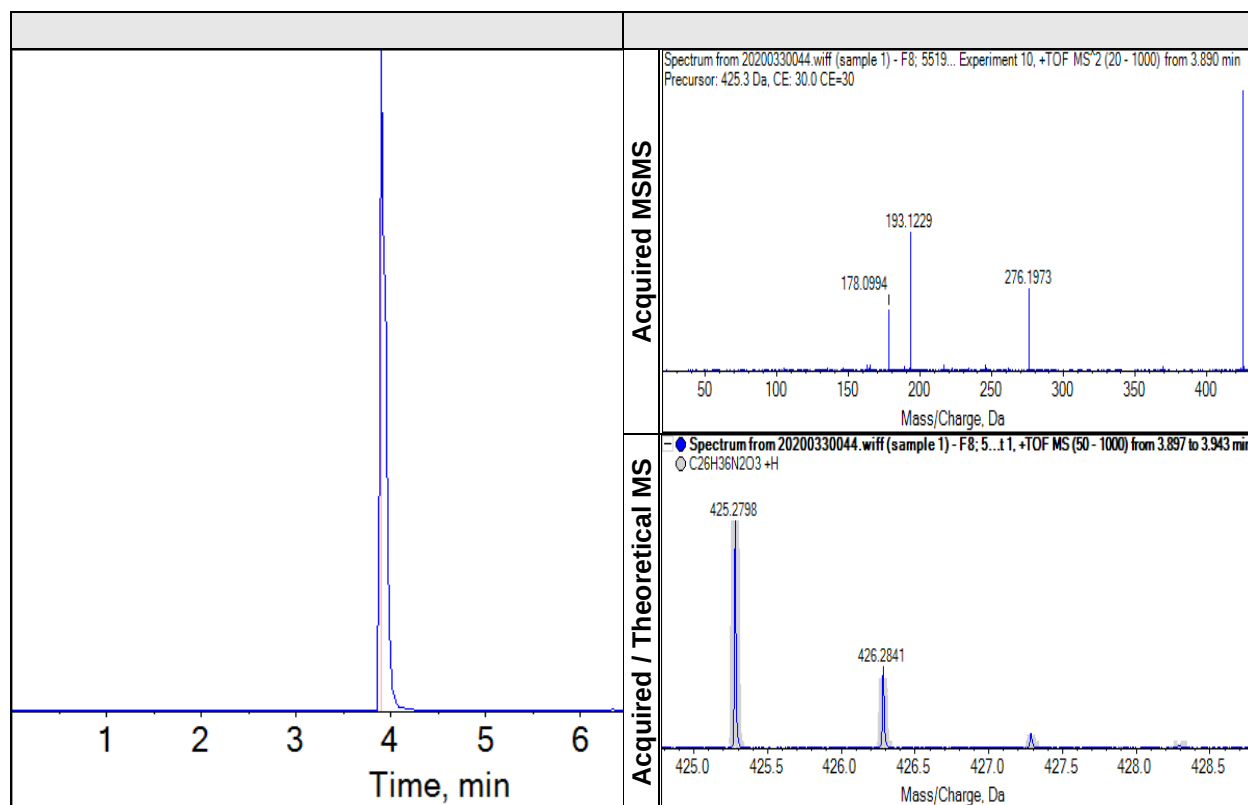


Figure 1: LC-QTOF Analysis of 28973

Table 1: LC-QTOF Conditions Used to Analyze 28973

Mass Spectrometer	Sciex TripleTOF 5600											
UPLC Hardware	Waters Acquity I-Class											
Mass Spectrometer Source	Turbolonspray, Positive											
Scan Ranges	TOF-MS: 50-1000 m/z MS/MS: 20-1000 m/z											
UPLC Column	Waters Acquity BEH C18, 2.1 × 50 mm, 1.7 μm											
Column Temperature	40°C											
Mobile Phase A	0.1% Formic Acid in Water											
Mobile Phase B	0.1% Formic Acid in Acetonitrile											
Gradient Table	<table><tr><td>Time, min</td><td>%B</td></tr><tr><td>0 – 0.5</td><td>5</td></tr><tr><td>0.5 – 5</td><td>Ramp 5 to 100</td></tr><tr><td>5 – 5.5</td><td>100</td></tr><tr><td>5.51 – 10</td><td>5</td></tr></table>		Time, min	%B	0 – 0.5	5	0.5 – 5	Ramp 5 to 100	5 – 5.5	100	5.51 – 10	5
Time, min	%B											
0 – 0.5	5											
0.5 – 5	Ramp 5 to 100											
5 – 5.5	100											
5.51 – 10	5											
Flow Rate	0.3 mL/min											
Injection Volume	5 μL											
Run Time	10 minutes											

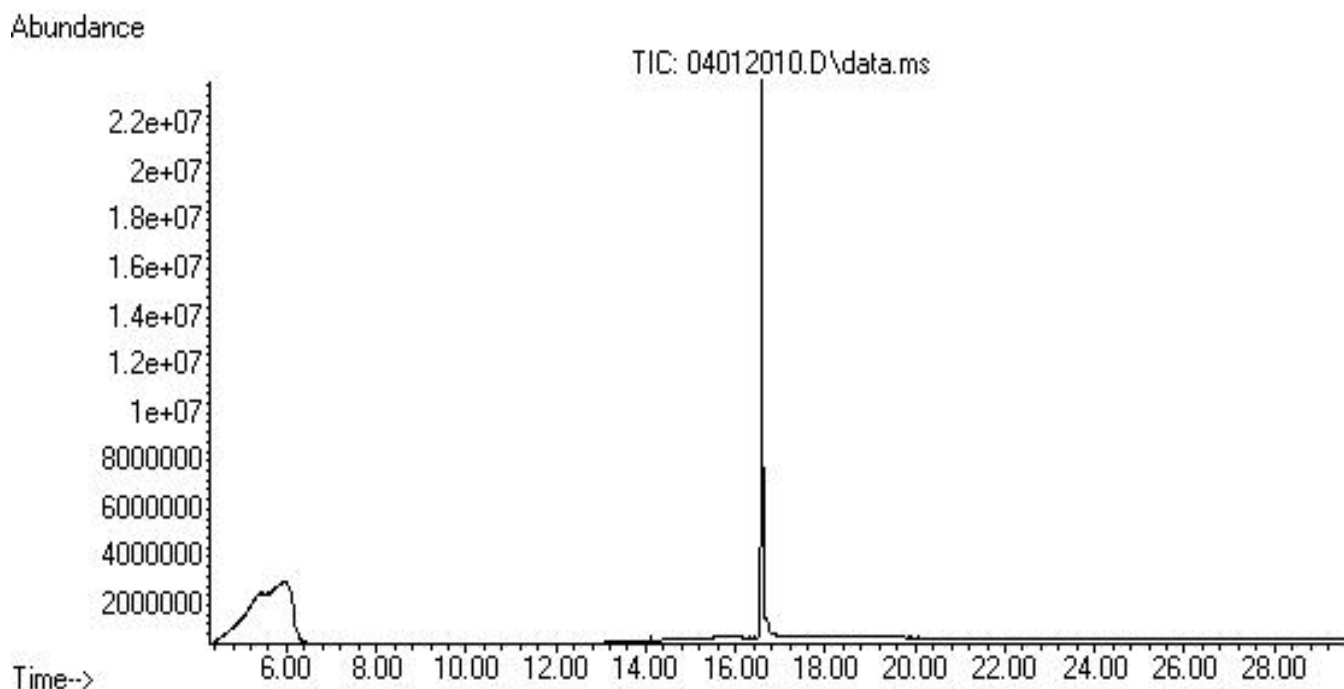
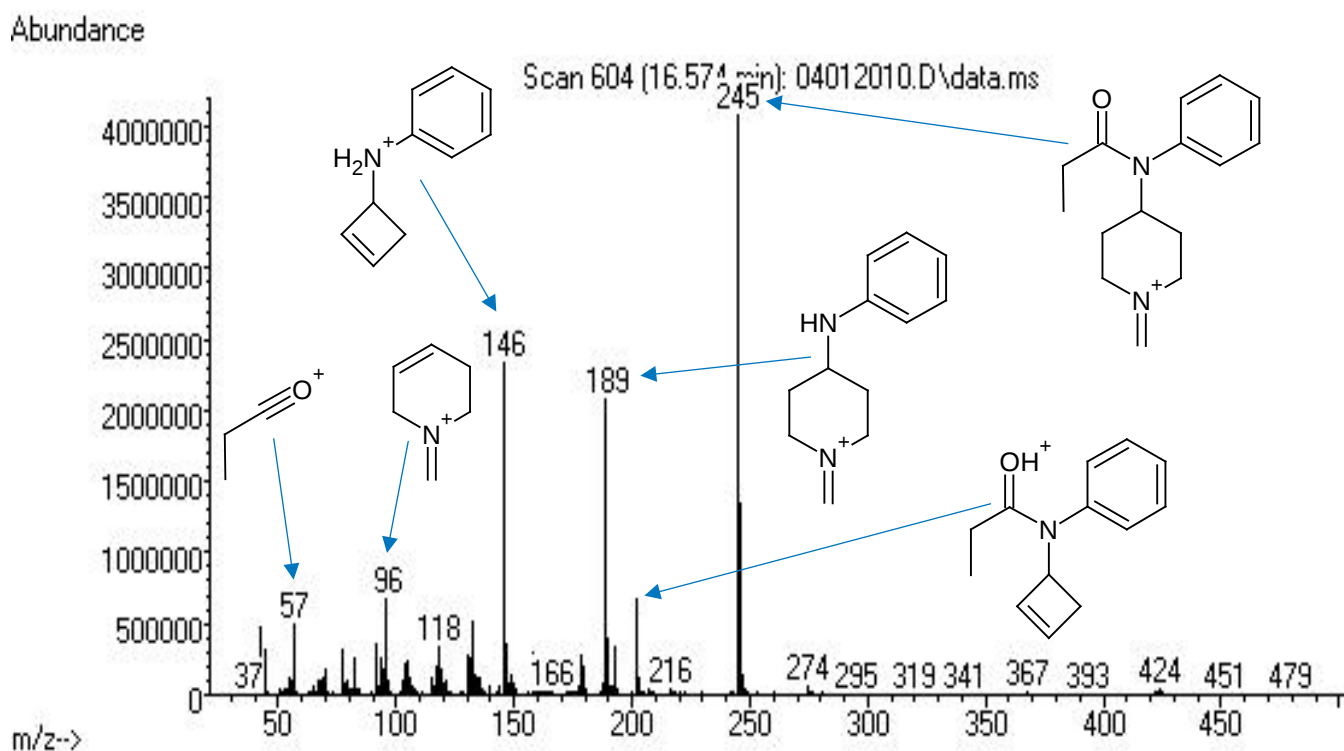


Figure 2. GC-MS Chromatogram (EI, TIC) of 28973



**Figure 3. Mass Spectrum (EI) of 28973 (Main Analyte)
Fragment Structure Assignments are Tentative and Unconfirmed**

Table 2. GC-MS Conditions Used to Analyze 28973

Instrument	Agilent 6890 GC/5973 MSD
Carrier Gas:	Helium
Column:	Restek Rtx-5MS 30m x 0.25mm x 0.25um
Oven Program:	45°C hold for 1.0 min 20°C/min. to 320°C, hold for 15 min.
Total Run Time:	30 minutes
Transfer Line Temp.:	280°C
Injection:	Cool-on-column
Injection Port Temperature:	Tracking oven temperature
Carrier Flow Rate:	1.6 mL/min.
Injection Volume:	0.5 uL
Acquisition Mode:	EI
Scan Range:	35 to 500 amu
Electron Impact:	70 eV
MS Quad Temperature:	150°C
MS Source Temperature:	230°C

F8; 55191-27-11

XEVO-TQXS#WBA0164

27-Mar-2020

27Mar20008 215 (5.833) Sm (Mn, 2x0.50)

3: Auto Daughters 425.26ES+

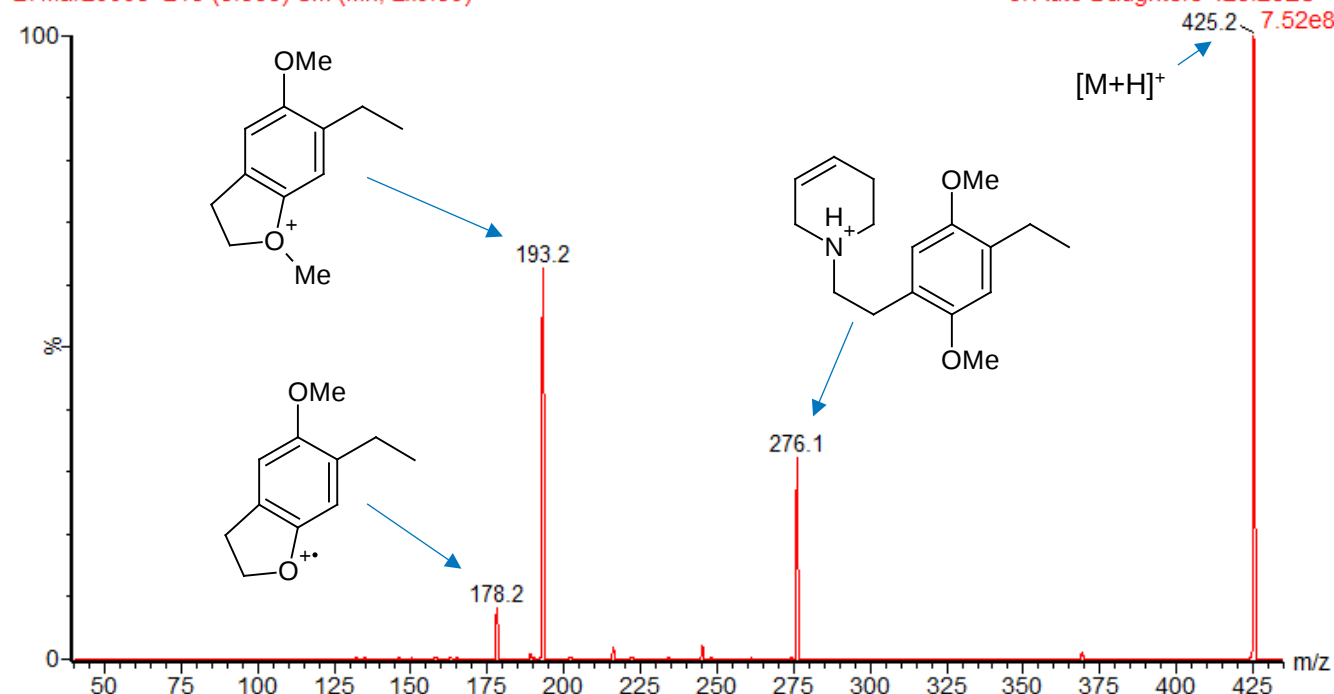


Figure 5. Low-resolution MS/MS (ESI, product ions of m/z 425) of 28973
Fragment Structure Assignments are Tentative and Unconfirmed

Table 3. LC-MS/MS-UV Conditions Used to Analyze 28973

Mass Spectrometer	Waters TQ-XS																	
UPLC Hardware	Waters Acquity I-Class																	
Mass Spectrometer Source	Electrospray, Positive Ionization																	
Photo Diode Array	Waters Acquity PDA																	
Scan Ranges	MS: 40-700 m/z PDA: 210-300 nm																	
Mass Selected for MS/MS fragmentation	Observed mass ± 0.25 Da																	
UPLC Column	Phenomenex Prodigy ODS (3), 100mm x 2.0mm, 3µm																	
Column Temperature	50°C																	
Mobile Phase A	1% Formic Acid in Water																	
Mobile Phase B	1% Formic Acid in Acetonitrile																	
Gradient Table	<table><tr><th>Time, min</th><th>%B</th><th>Flow, mL/min</th></tr><tr><td>0 – 1</td><td>2</td><td>0.3</td></tr><tr><td>1 – 11</td><td>Ramp 2 to 100</td><td>0.3</td></tr><tr><td>11.1 – 12</td><td>100</td><td>0.5</td></tr><tr><td>12.1 – 14</td><td>2</td><td>0.3</td></tr></table>			Time, min	%B	Flow, mL/min	0 – 1	2	0.3	1 – 11	Ramp 2 to 100	0.3	11.1 – 12	100	0.5	12.1 – 14	2	0.3
Time, min	%B	Flow, mL/min																
0 – 1	2	0.3																
1 – 11	Ramp 2 to 100	0.3																
11.1 – 12	100	0.5																
12.1 – 14	2	0.3																
Injection Volume	25 µL																	
Run Time	14 minutes																	