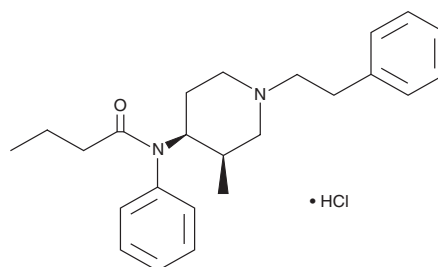


# CERTIFICATE of ANALYSIS

(±)-*cis*-3-methyl Butyryl fentanyl  
(hydrochloride)



ACCREDITED  
ISO/IEC 17025 #AT-1773  
ISO 17034 #AR-1774

|                       |                                                                                                                                                                                                                                                                                                                                                                                   |
|-----------------------|-----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Item No.:             | 20817                                                                                                                                                                                                                                                                                                                                                                             |
| Batch No.:            | 0560036                                                                                                                                                                                                                                                                                                                                                                           |
| CAS Registry No.:     | 88641-20-7                                                                                                                                                                                                                                                                                                                                                                        |
| Molecular Formula:    | C <sub>24</sub> H <sub>32</sub> N <sub>2</sub> O • HCl                                                                                                                                                                                                                                                                                                                            |
| Formula Weight:       | 401.0 amu                                                                                                                                                                                                                                                                                                                                                                         |
| UV λ <sub>max</sub> : | 208, 258 nm                                                                                                                                                                                                                                                                                                                                                                       |
| Expiry Date:          | 16JUL2026 (valid from date of certification)                                                                                                                                                                                                                                                                                                                                      |
| Supplied as:          | A neat solid                                                                                                                                                                                                                                                                                                                                                                      |
| Storage:              | Unopened at -20°C                                                                                                                                                                                                                                                                                                                                                                 |
| Safety:               | Refer to Safety Data Sheet                                                                                                                                                                                                                                                                                                                                                        |
| Intended Use:         | For analytical testing purposes only, not intended for human or animal use.                                                                                                                                                                                                                                                                                                       |
| Instructions for Use: | Once opened this material should be minimally exposed to ambient conditions and returned to recommended storage conditions immediately after use. Ongoing stability testing supports a negligible decrease in purity over a series of thaw-refreeze cycles. It is recommended that laboratories perform periodic testing to verify the material remains fit for the intended use. |

Approval: 

Title: ISO Quality Manager

Certification Date: 16JUL2019

Cayman Chemical certifies that this standard meets the specifications stated in this certificate and warrants this product to meet the stated acceptance criteria through the expiration date when stored unopened as recommended.



# CERTIFICATE of ANALYSIS

| Qualifier                    | Method                       | Result              |
|------------------------------|------------------------------|---------------------|
| Appearance                   | Visual inspection            | White solid         |
| Chromatographic Purity, HPLC | Cayman Method TST SD132      | >99.90% $\pm$ 0.18% |
| Identity, LC-MS              | Cayman Method TST SD13, +ESI | 365.2 amu           |
| Identity, GC-MS              | Cayman Method TST SD12       | Conforms            |
| Identity, FTIR               | USP<854> (TST SD03)          | Conforms            |
| Identity, NMR                | $^1\text{H}$ NMR             | Conforms            |

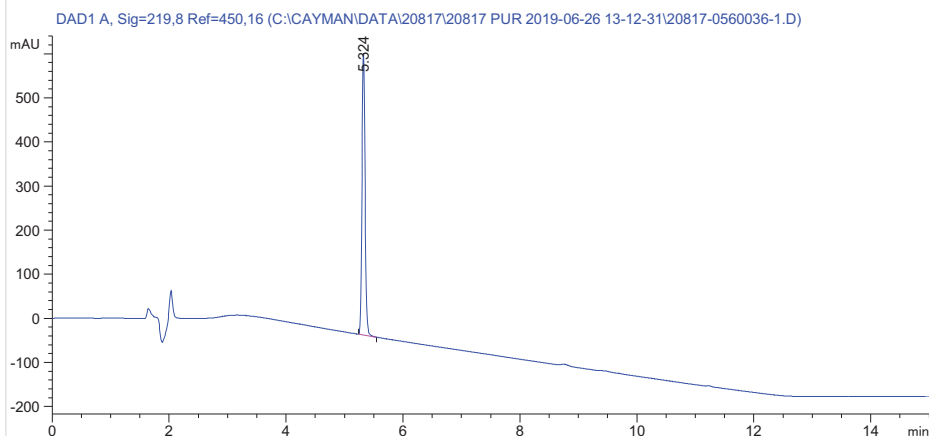
NMR and optical rotation (if applicable) are provided as supplemental information but are not within scope of ISO accreditation. Property values are traceable to SI units through an unbroken chain of measurements.

## Measurement Uncertainty

All measurement uncertainties are expressed as expanded uncertainties in accordance with ISO standards for Testing Laboratories and Reference Material Producers at the approximate 95% confidence interval using an appropriate coverage factor. Where applicable, optical rotation, chiral purity, and/or isotopic purity testing are performed to support the identification of the reference material, therefore the uncertainty is considered null.

## Supplemental Data (Neat Material)

### HPLC-UV

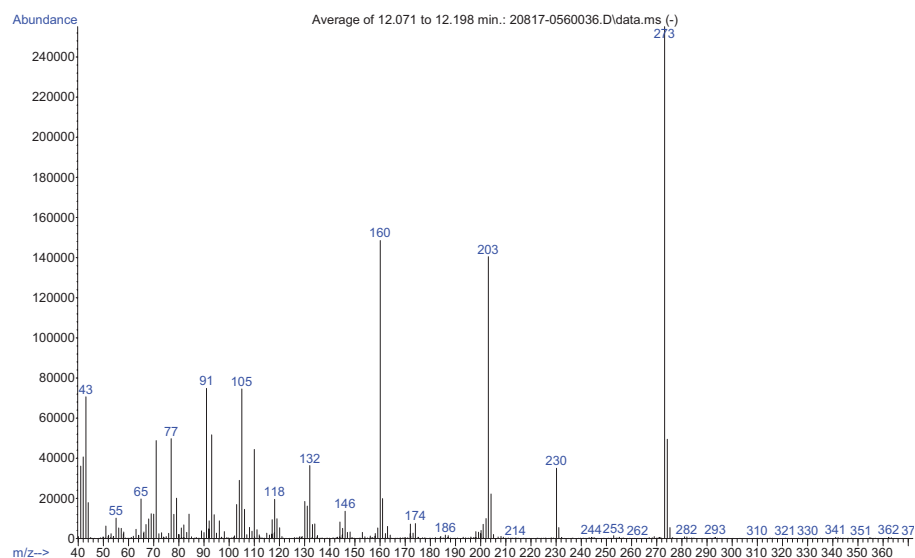


### Conditions

| Instrument   | Agilent 1100/1200 Series                                                                                                                                                                                                          |            |    |      |        |       |     |         |     |
|--------------|-----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|------------|----|------|--------|-------|-----|---------|-----|
| Column       | 4.6 x 150 mm, 5 $\mu\text{m}$ Zorbax                                                                                                                                                                                              |            |    |      |        |       |     |         |     |
| Mobile Phase | A: 0.1% Trifluoroacetate<br>B: Acetonitrile                                                                                                                                                                                       |            |    |      |        |       |     |         |     |
| Gradient     | <table border="1"> <thead> <tr> <th>Time (min)</th> <th>%B</th> </tr> </thead> <tbody> <tr> <td>0-10</td> <td>20-95%</td> </tr> <tr> <td>10-13</td> <td>95%</td> </tr> <tr> <td>13.1-20</td> <td>20%</td> </tr> </tbody> </table> | Time (min) | %B | 0-10 | 20-95% | 10-13 | 95% | 13.1-20 | 20% |
| Time (min)   | %B                                                                                                                                                                                                                                |            |    |      |        |       |     |         |     |
| 0-10         | 20-95%                                                                                                                                                                                                                            |            |    |      |        |       |     |         |     |
| 10-13        | 95%                                                                                                                                                                                                                               |            |    |      |        |       |     |         |     |
| 13.1-20      | 20%                                                                                                                                                                                                                               |            |    |      |        |       |     |         |     |
| Flow Rate    | 1.0 ml/min                                                                                                                                                                                                                        |            |    |      |        |       |     |         |     |
| Column Temp  | 30°C                                                                                                                                                                                                                              |            |    |      |        |       |     |         |     |
| Wavelength   | UV monitored at 219 nm                                                                                                                                                                                                            |            |    |      |        |       |     |         |     |

# CERTIFICATE of ANALYSIS

## GC-MS

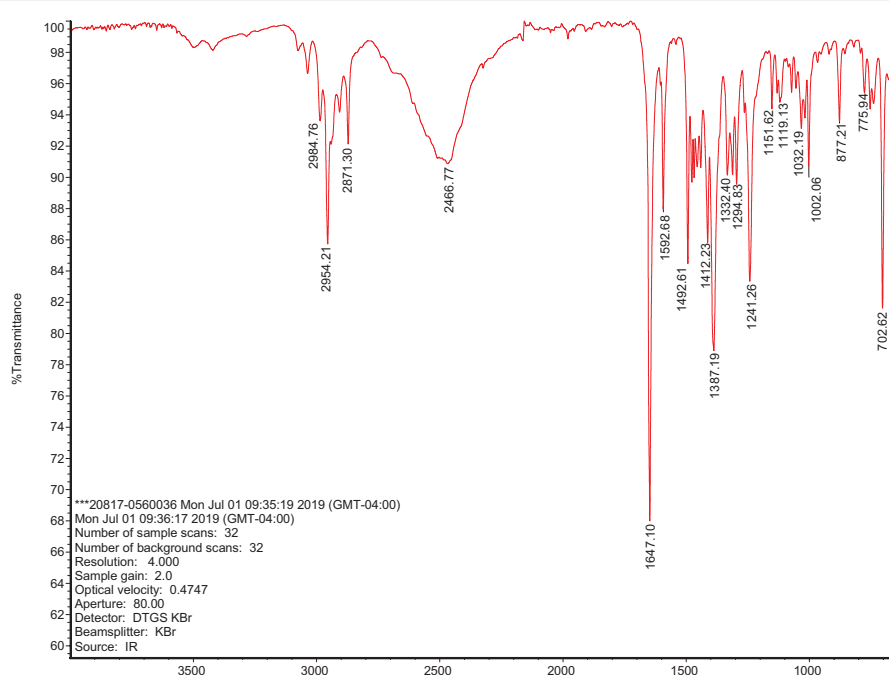


## Conditions

|                    |                                                                                   |
|--------------------|-----------------------------------------------------------------------------------|
| Instrument         | Agilent 6890 GC/5973 MSD                                                          |
| Column             | 30 m x 0.32 mm,<br>0.5 µm Rtx-5MS                                                 |
| Carrier Gas        | He                                                                                |
| Flow Rate          | 2 ml/min                                                                          |
| Inlet Temp         | 300°C                                                                             |
| Split Ratio        | 15:1                                                                              |
| Oven Program       | 50°C hold for 1 min, ramp<br>to 300°C at 30°C per min,<br>hold at 300°C to 15 min |
| Transfer Line Temp | 300°C                                                                             |
| Voltage            | 70eV EI MS                                                                        |
| Scan Range         | 40-600 m/z                                                                        |
| Tune File          | stune                                                                             |

Apex spectrum – background (1 min window in front of peak)

## FTIR



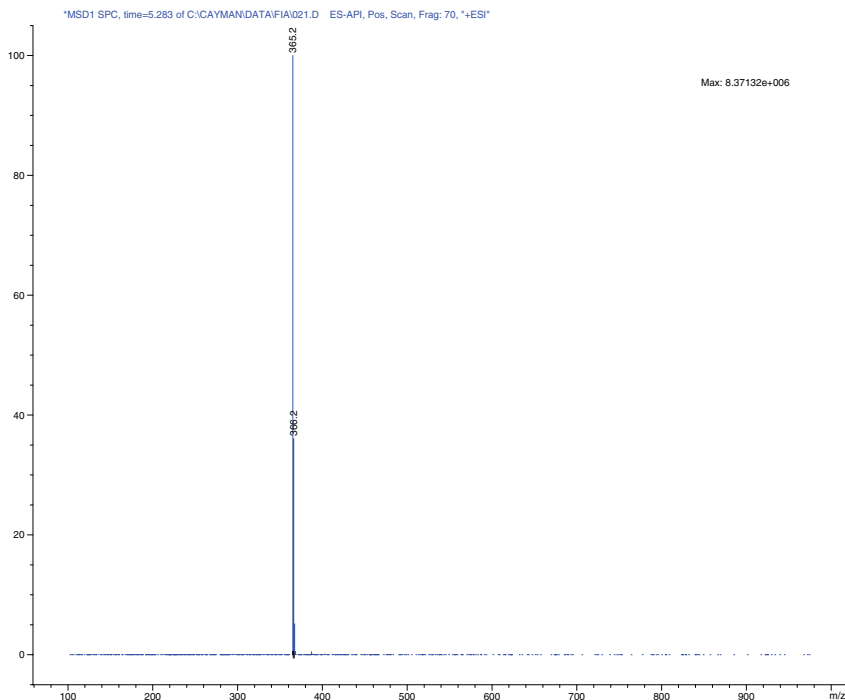
## Conditions

|            |                                                                   |
|------------|-------------------------------------------------------------------|
| Instrument | Thermo Nicolet iS10 FTIR /<br>Diamond SmartATR (single<br>bounce) |
| Scans      | 32 scans /<br>32 background scans                                 |
| Range      | 650-4,000 cm <sup>-1</sup>                                        |
| Resolution | 4.000                                                             |

ATR and background corrected

# CERTIFICATE of ANALYSIS

## ESI-MS



## Conditions

|                      |                                                |
|----------------------|------------------------------------------------|
| Instrument           | Agilent HPLC/MSD                               |
| Mobile Phase         | 50:50:0.1<br>MeOH/H <sub>2</sub> O/Acetic acid |
| Flow Rate            | 0.5 ml/min                                     |
| Ionization Mode      | +ESI                                           |
| Mass Range           | 100-1,000 m/z                                  |
| Nebulizer            | 60 psi                                         |
| Desolvation Gas      | 13 L/min                                       |
| Desolvation Temp     | 350°C                                          |
| Electrospray Voltage | 4kV                                            |

MS collected across peak width at half height

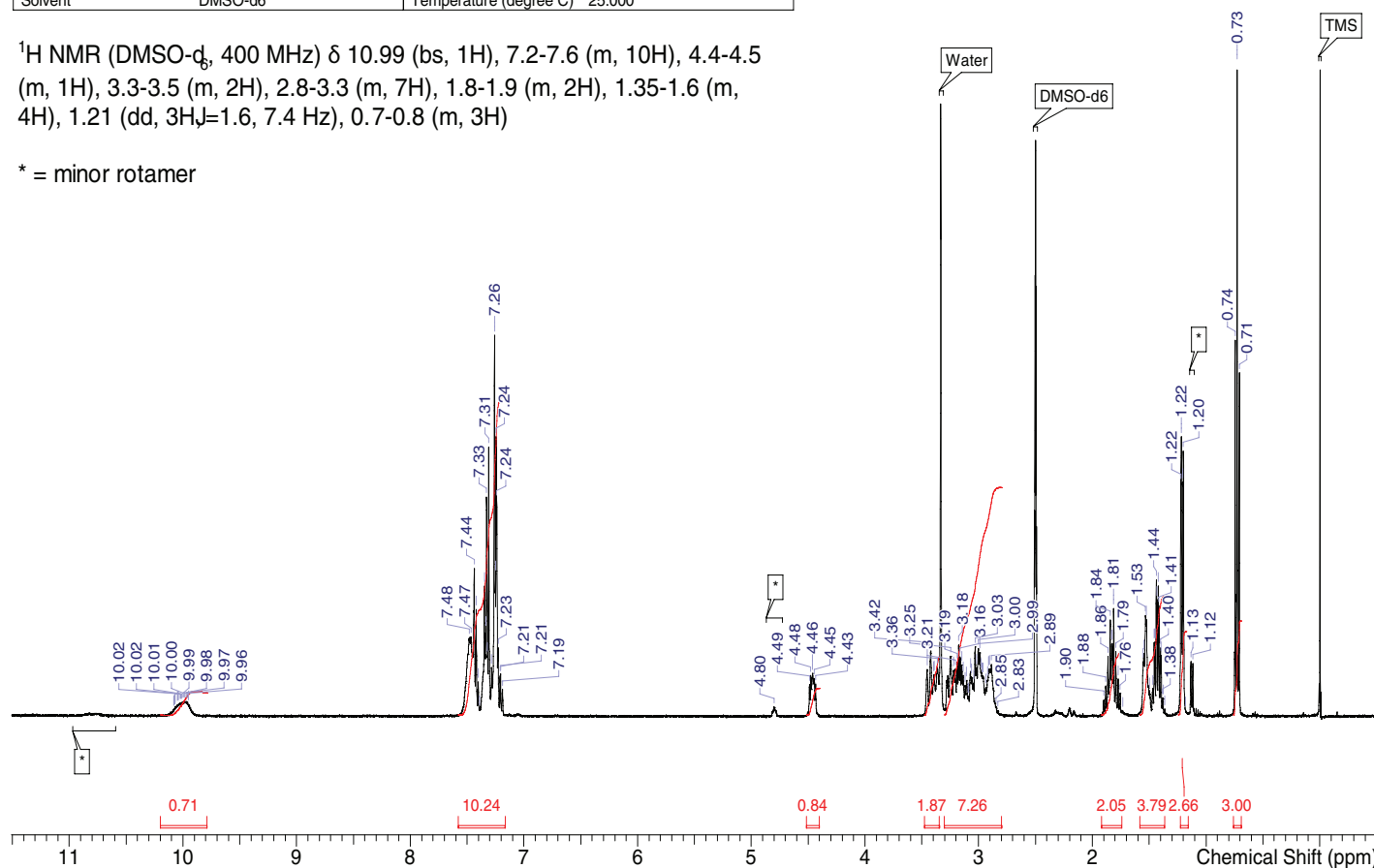
# CERTIFICATE of ANALYSIS

## NMR (not within scope of ISO accreditation)

|                 |                                                                                                                      |                        |        |                         |
|-----------------|----------------------------------------------------------------------------------------------------------------------|------------------------|--------|-------------------------|
| Date            | Jun 24 2019                                                                                                          |                        |        |                         |
| File Name       | \\sulfur\private\nmrdata\ 2019\20817-0560036_20190624_1\20817-0560036_20190624_01\20817-0560036_20190624_001.fid\fid |                        |        |                         |
| Frequency (MHz) | 399.9680                                                                                                             | Nucleus                | 1H     | Number of Transients 32 |
| Solvent         | DMSO-d6                                                                                                              | Temperature (degree C) | 25.000 |                         |

<sup>1</sup>H NMR (DMSO-d<sub>6</sub>, 400 MHz) δ 10.99 (bs, 1H), 7.2-7.6 (m, 10H), 4.4-4.5 (m, 1H), 3.3-3.5 (m, 2H), 2.8-3.3 (m, 7H), 1.8-1.9 (m, 2H), 1.35-1.6 (m, 4H), 1.21 (dd, 3H, J=1.6, 7.4 Hz), 0.7-0.8 (m, 3H)

\* = minor rotamer



### Conditions

|            |                          |
|------------|--------------------------|
| Instrument | Varian Inova 400 MHz NMR |
| Scans      | 64 scans                 |

### Homogeneity

A minimum sample size of 2.0 µg was used to determine homogeneity of the bulk solid. The recommended minimum quantity for use is 2.0 µg. Quantities below this have not been evaluated.

### Short-Term Stability

No decrease in the purity was observed at ambient or 60°C after two weeks. This data supports shipping of this product at ambient temperature.

### Long-Term Stability

Long-term stability data predicts seven years stability at the -20°C storage temperature. Long-term stability studies are ongoing and the Certificate of Analysis will be updated upon study completion.



# CERTIFICATE of ANALYSIS

## Revision History

| Revision No. | Date      | Reason for Revision |
|--------------|-----------|---------------------|
| 01           | 16JUL2019 | Initial version     |

## Disclaimers

### Material Safety Data

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some but not all of the information required for the safe and proper use of this material. Before use, review the complete Safety Data Sheet, which has been sent *via* email to your institution.

### Warranty and Limitation of Remedy

Cayman Chemical Company makes no warranty or guarantee of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specification.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver of Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitations of Remedy located on our website and in our catalog.

This Certificate shall not be reproduced except in full, without written approval from the Cayman Chemical ISO Quality Manager.

ISO CRT SD01 v 3.1

**CAYMAN CHEMICAL**  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335  
**FAX:** [734] 971-3640

crmquality@caymanchem.com  
www.caymanchem.com