

CERTIFICATE OF ANALYSIS

4'-fluoro, para-fluoro (±)-trans-3-methyl Fentanyl (hydrochloride)

N-((3S,4S)-1-(4-fluorophenethyl)-3-methylpiperidin-4-yl)-N-(4-fluorophenyl)propionamide, monohydrochloride

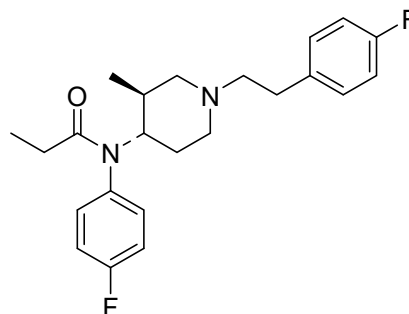
Item #: 24503
Batch #: 0528336
CAS Registry Number: N/A

Expiry Date: 24APR2023 (valid from date of certification)

Description: neat solid

Storage and Handling: Store at -20°C. Warm to room temperature prior to opening.

Safety: Poison



Compound Information

Chemical Formula	C ₂₃ H ₂₈ F ₂ N ₂ O • HCl
Formula Weight	422.9 amu
UV λ _{max}	210 nm



ISO/IEC 17025
#AT-1773
ISO Guide 34
#AR-1774



Quality Information

Qualifier	Method	Limit	Result	Meets Specification
Appearance	Visual inspection	White/off-white solid	White solid	Y
Chromatographic Purity, HPLC	Cayman Method TST SD132	≥98.00%	99.41% ± 0.18%	Y
Identity, LCMS	Cayman Method TST SD13, +ESI	387.2 ± 0.5 amu	387.2 amu	Y
Identity, GCMS	Cayman Method TST SD12	Conforms	Conforms	Y
FTIR	USP<851> (diamond ATR)	Conforms	Conforms	Y
*Identity, NMR	¹ H NMR	Conforms	Conforms	Y

*NMR is provided as supplemental info but is 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 17025 and Guide 34 at the approximate 95% confidence interval using the appropriate coverage factor.

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.

Approval:  Title: Cayman Chemical ISO Quality Manager

Certification Date: 24APR2018

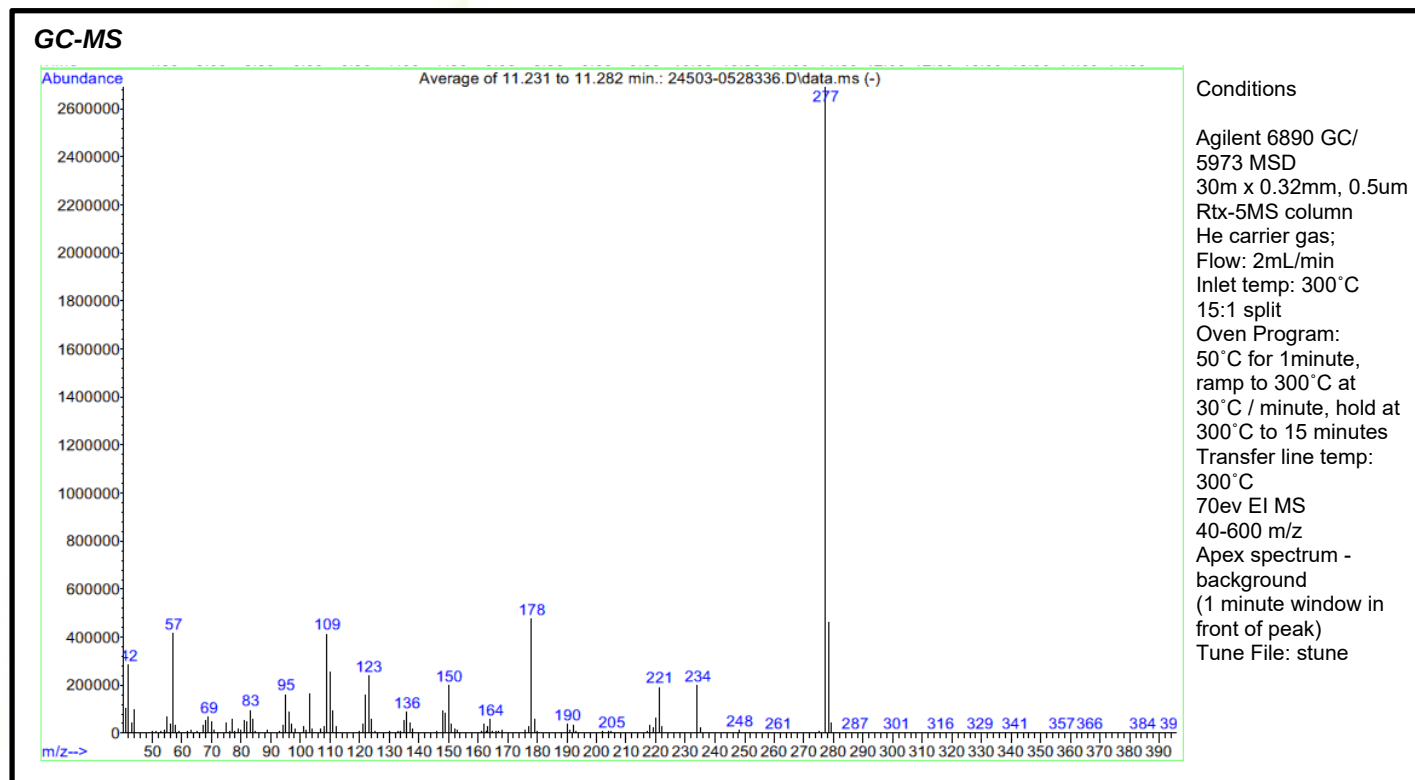
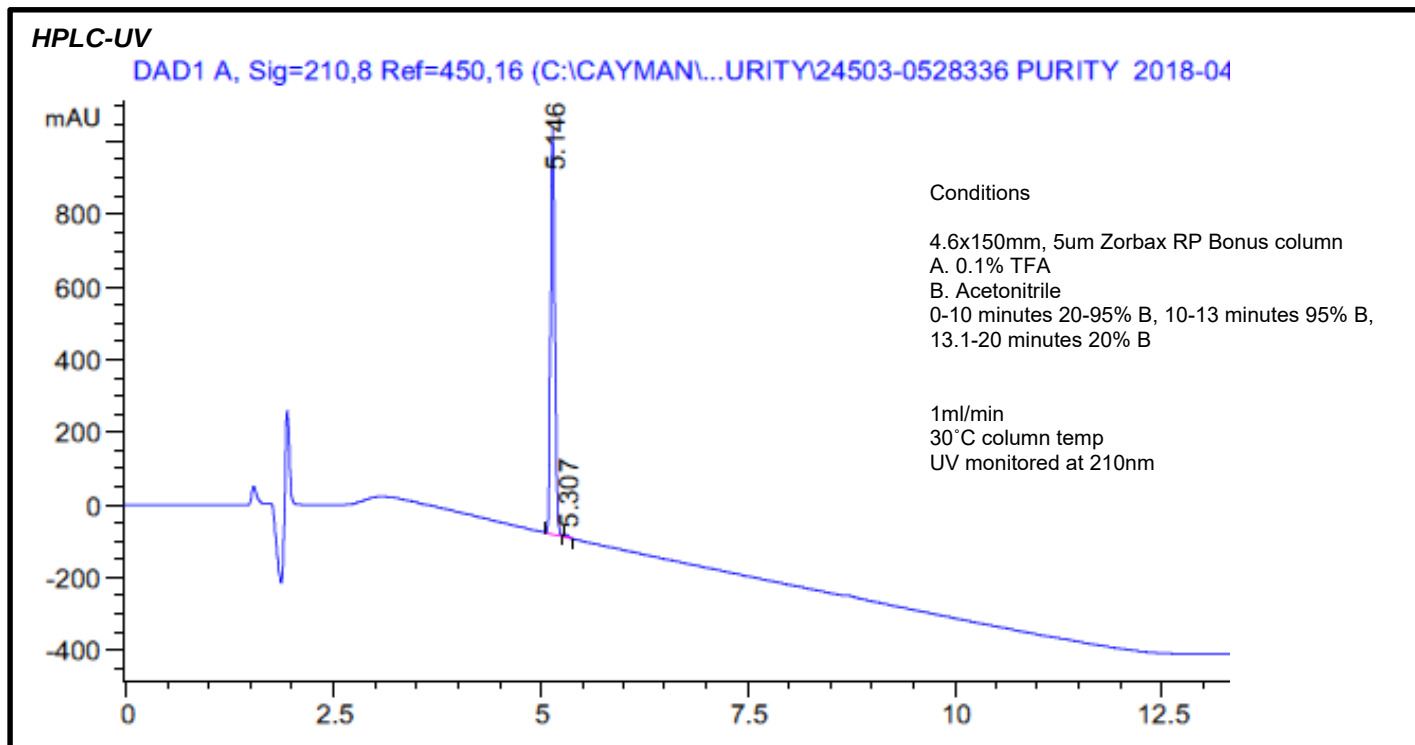
Cayman Chemical Co.

1180 E. Ellsworth Rd
Ann Arbor, MI 48108
www.caymanchem.com

Contact Us:

1-800-364-9897 or (734)971-3335
Fax: (734)971-3640
Email: crmquality@caymanchem.com

For R&D purposes only. Not intended for human or animal use.

Supplemental Data

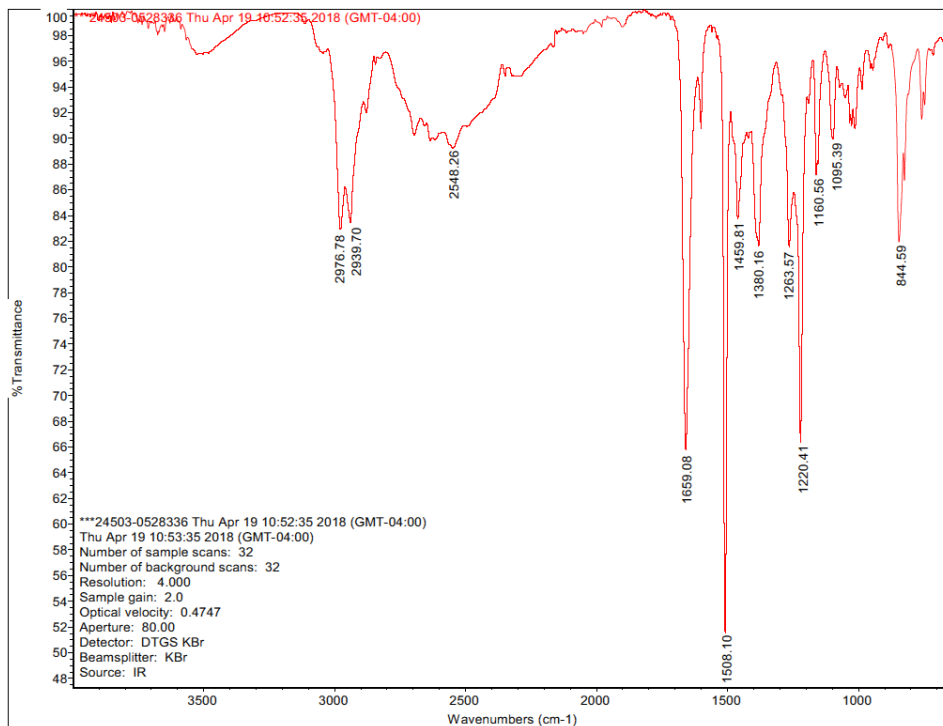
Cayman Chemical Co.

1180 E. Ellsworth Rd
Ann Arbor, MI 48108
www.caymanchem.com

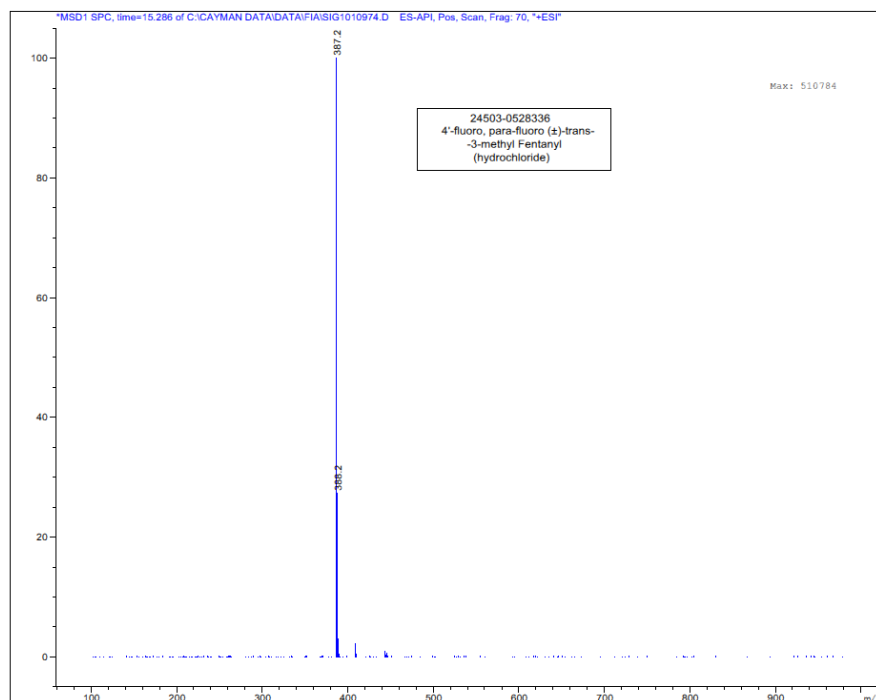
Contact Us:

1-800-364-9897 or (734)971-3335
Fax: (734)971-3640
Email: crmquality@caymanchem.com

For R&D purposes only. Not intended for human or animal use.

FTIR**Conditions**

Thermo Nicolet iS10 FTIR /
Diamond SmartATR (single bounce)
32 scans / 32 background scans
Range: 650-4000cm⁻¹
Resolution: 4.000
ATR and background corrected

ESI-MS**Conditions**

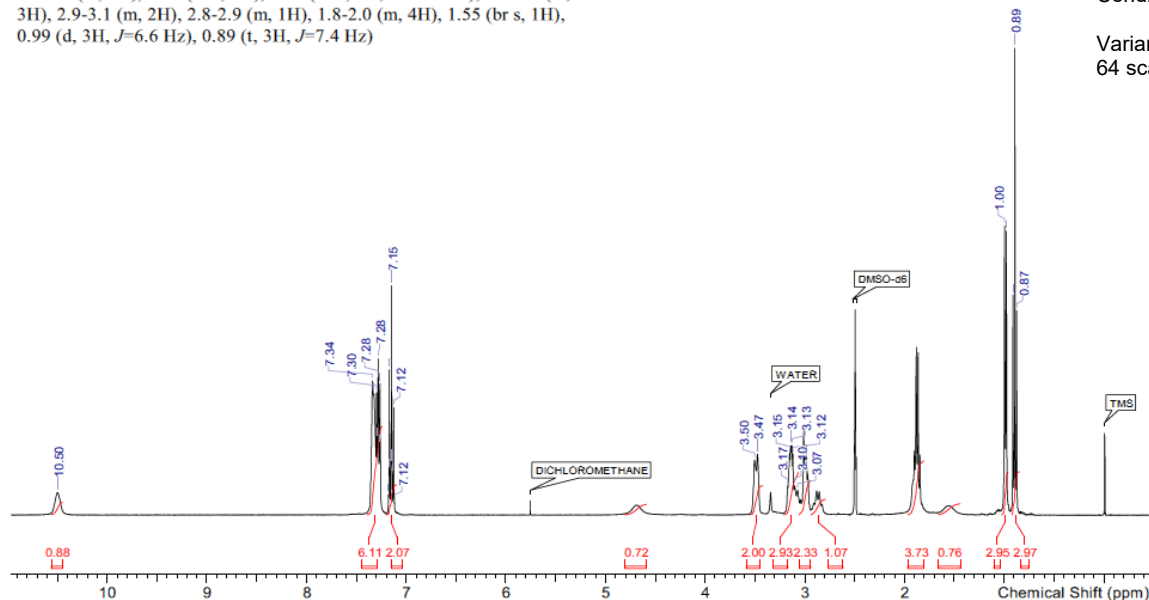
Agilent 1100 HPLC / G1956MSD
FIA-MS (no column)
Mobile phase: 50:50:0.1
MeOH/H₂O/acetic acid
Flow: 0.5mL/min
+ESI 100-1000m/z range
Nebulizer: 60psi
Desolvation Gas: 13L/min
Desolvation Temp: 350°C
Electrospray Voltage: 4kV
MS collected across peak width
at half height

NMR *not within scope of ISO 17025/Guide 34 accreditation

Date	Apr 5 2018
File Name	\\sulfur\private\nmrdata\ 2018\24503-0528336_20180405_124503-0528336_20180405_01\24503-0528336_20180405_001.fid\fid
Frequency (MHz)	399.9680
Solvent	DMSO-d6
Nucleus	1H
Temperature (degree C)	25.000
Number of Transients	32

¹H NMR (DMSO-d₆, 400 MHz) δ 10.50 (br s, 1H), 7.2-7.4 (m, 6H), 7.1-7.2 (m, 2H), 4.70 (br s, 1H), 3.49 (br d, 2H, *J*=11.3 Hz), 3.1-3.2 (m, 3H), 2.9-3.1 (m, 2H), 2.8-2.9 (m, 1H), 1.8-2.0 (m, 4H), 1.55 (br s, 1H), 0.99 (d, 3H, *J*=6.6 Hz), 0.89 (t, 3H, *J*=7.4 Hz)

Conditions

Varian Inova 400MHz NMR
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 Summary

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

Long Term Stability

Long term stability data predicts 5 years stability at the -20°C storage temperature.

Cayman Chemical Co.

1180 E. Ellsworth Rd
Ann Arbor, MI 48108
www.caymanchem.com

Contact Us:

1-800-364-9897 or (734)971-3335
Fax: (734)971-3640
Email: crmquality@caymanchem.com

For R&D purposes only. Not intended for human or animal use.

Revision History

Revision #	Date	Reason for Revision
01	24APR2018	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 Material 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 1.1

