

# CERTIFICATE of ANALYSIS



## para-methyl Furanyl fentanyl (hydrochloride)

N-(1-phenethylpiperidin-4-yl)-N-(p-tolyl)furan-2-carboxamide, monohydrochloride

Item No. 22713 • Batch No. 0511784

Purity Specification:  $\geq 98\%$

Molecular Formula : C<sub>25</sub>H<sub>28</sub>N<sub>2</sub>O<sub>2</sub> • HCl

CAS Number: (unspecified)

Formula Weight : 425

Expiry date: 4/30/2021

| Tests         | Results               |
|---------------|-----------------------|
| HPLC          | Purity: 100.0 %       |
| IR            | Conforms              |
| Mass spec     | M-HCl: 389.2          |
| Melting Point | 260 - 262 °C          |
| TLC           | Purity: 100 %         |
| UV            | $\lambda$ max: 258 nm |
| NMR           | Conforms              |

Reviewed and approved by: Jennifer LaBrecque

A handwritten signature in black ink that reads "Jen LaBrecque".

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 • USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM