

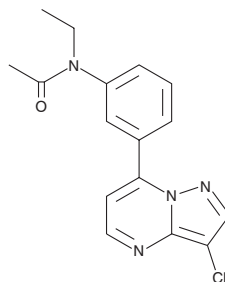
PRODUCT INFORMATION



Zaleplon

Item No. 11577

CAS Registry No.: 151319-34-5
Formal Name: N-[3-(3-cyanopyrazolo[1,5-a]pyrimidin-7-yl)phenyl]-N-ethyl-acetamide
Synonym: CL 284,846
MF: C₁₇H₁₅N₅O
FW: 305.3
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥1 year



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Zaleplon (Item No. 11577) is an analytical reference material that is functionally categorized as a sedative. It is a pyrazolopyrimidine that selectively activates the GABA_A α₁ receptor subunit (EC₅₀ = 1.1 μM), producing a sedative effect.¹⁻³ It is characterized as being an ultra-short-acting sedative due to short T_{max} (0.7-1.4 hours) and half-life (1 hour) values.² This product is intended for research and forensic applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards.

References

1. Petroski, R.E., Pomeroy, J.E., Das, R., *et al.* Indiplon is a high-affinity positive allosteric modulator with selectivity for α₁ subunit-containing GABA_A receptors. *J. Pharmacol. Exp. Ther.* **317**(1), 369-377 (2006).
2. Gunja, N. The clinical and forensic toxicology of Z-drugs. *J. Med. Toxicol.* **9**, 155-162 (2013).
3. Nutt, D.J., Stahl, S.M. Searching for perfect sleep: The continuing evolution of GABA_A receptor modulators as hypnotics. *J. Psychopharmacol.* **24**(11), 1601-1612 (2010).

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, 08/08/2019

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