

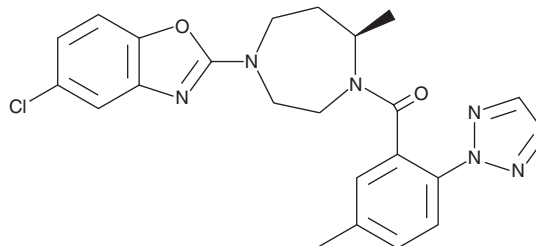
PRODUCT INFORMATION



Suvorexant

Item No. 22911

CAS Registry No.: 1030377-33-3
Formal Name: [(7R)-4-(5-chloro-2-benzoxazolyl)hexahydro-7-methyl-1H-1,4-diazepin-1-yl][5-methyl-2-(2H-1,2,3-triazol-2-yl)phenyl]-methanone
Synonym: MK-4305
MF: C₂₃H₂₃ClN₆O₂
FW: 450.9
Purity: ≥98%
UV/Vis.: λ_{max}: 216, 258 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

Suvorexant (Item No. 22911) is an analytical reference standard categorized as a diazepane sedative.¹ Formulations containing suvorexant show abuse potential similar to zolpidem (Item No. 15792) in recreational polydrug users with a history of sedative and psychedelic drug use.² Suvorexant is regulated as a Schedule IV compound in the United States. This product is intended for use in analytical forensic applications.

References

1. Winrow, C.J. and Renger, J.J. Discovery and development of orexin receptor antagonists as therapeutics for insomnia. *Br. J. Pharmacol.* **171**(2), 283-293 (2014).
2. Schoedel, K.A., Sun, H., Sellers, E.M., *et al.* Assessment of the abuse potential of the orexin receptor antagonist, suvorexant, compared with zolpidem in a randomized crossover study. *J. Clin. Psychopharmacol.* **36**(4), 314-323 (2016).

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, 06/16/2023

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