

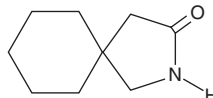
# PRODUCT INFORMATION



## Gabapentin lactam

Item No. 43246

CAS Registry No.: 64744-50-9  
Formal Name: 2-azaspiro[4.5]decan-3-one  
MF:  $C_9H_{15}NO$   
FW: 153.2  
Purity:  $\geq 98\%$   
Supplied as: A neat solid  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 4$  years



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

### Description

Gabapentin lactam (Item No. 43246) is an analytical reference standard categorized as a degradation product of gabapentin.<sup>1,2</sup> Gabapentin lactam is a potential impurity in gabapentin preparations.<sup>1</sup> It has convulsant activity in rats.<sup>3</sup> This product is intended for research and forensic applications.

### References

1. Ciavarella, A.B., Gupta, A., Sayeed, V.A., *et al.* Development and application of a validated HPLC method for the determination of gabapentin and its major degradation impurity in drug products. *J. Pharm. Biomed. Anal.* **43**(5), 1647-1653 (2007).
2. He, Y., Jia, D., Du, S., *et al.* Toxicity of gabapentin-lactam on the early developmental stage of zebrafish (*Danio rerio*). *Environ. Pollut.* **287**, 117649 (2021).
3. Potschka, H., Feuerstein, T.J., and Löscher, W. Gabapentin-lactam, a close analogue of the anticonvulsant gabapentin, exerts convulsant activity in amygdala kindled rats. *Naunyn Schmiedebergs Arch. Pharmacol.* **361**(2), 200-205 (2000).

#### 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, 04/24/2025

#### 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