

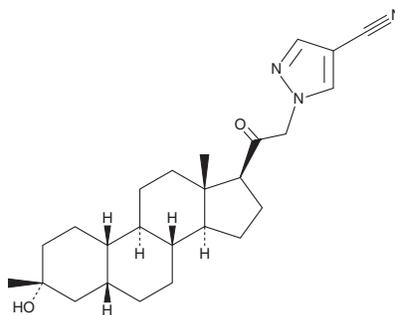
# PRODUCT INFORMATION



## SAGE-217

Item No. 31755

**CAS Registry No.:** 1632051-40-1  
**Formal Name:** 1-[(3 $\alpha$ ,5 $\beta$ )-3-hydroxy-3-methyl-20-oxo-19-norpregnan-21-yl]-1H-pyrazole-4-carbonitrile  
**Synonyms:** S-812217, SGE 797, Zuranolone  
**MF:** C<sub>25</sub>H<sub>35</sub>N<sub>3</sub>O<sub>2</sub>  
**FW:** 409.6  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\max}$ : 215 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



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

### Description

SAGE-217 (Item No. 31755) is an analytical reference standard categorized as a neuroactive steroid and GABA<sub>A</sub> receptor positive allosteric modulator.<sup>1,2</sup> SAGE-217 has anticonvulsant activity in mice and reduces neuropathic pain in a mouse model of segmental spinal nerve ligation.<sup>2,3</sup> Formulations containing SAGE-217 have been used in the treatment of postpartum depression. SAGE-217 is regulated as a Schedule IV compound in the United States. This product is intended for research and forensic applications.

### References

1. Botella, G.M., Salituro, F.G., Harrison, B.L., *et al.* Neuroactive steroids. 2. 3 $\alpha$ -Hydroxy-3 $\beta$ -methyl-21-(4-cyano-1H-pyrazol-1'-yl)-19-nor-5 $\beta$ -pregnan-20-one (SAGE-217): A clinical next generation neuroactive steroid positive allosteric modulator of the ( $\gamma$ -aminobutyric acid)<sub>A</sub> receptor. *J. Med. Chem.* **60**(18), 7810-7819 (2017).
2. Althaus, A.L., Ackley, M.A., Belfort, G.M., *et al.* Preclinical characterization of zuranolone (SAGE-217), a selective neuroactive steroid GABA<sub>A</sub> receptor positive allosteric modulator. *Neuropharmacology* **181**, 108333 (2020).
3. Xu, C., Wu, J.H., Yu, H., *et al.* Effect of two novel GABAA receptor positive allosteric modulators on neuropathic and inflammatory pain in mice. *Neuropharmacology* **269**, 110317 (2025).

#### 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, 08/25/2025

#### 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