

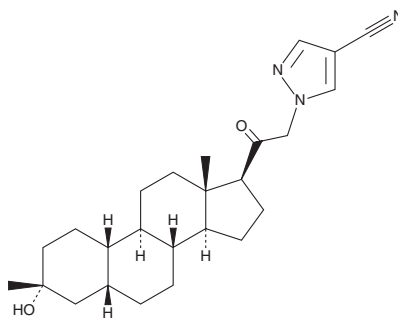
# PRODUCT INFORMATION



## 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.:**  $\lambda_{\text{max}}$ : 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.

### Laboratory Procedures

SAGE-217 is supplied as a crystalline solid. A stock solution may be made by dissolving the SAGE-217 in the solvent of choice, which should be purged with an inert gas. SAGE-217 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of SAGE-217 is approximately 10 mg/ml in DMSO and 30 mg/ml in DMF.

SAGE-217 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SAGE-217 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. SAGE-217 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

SAGE-217 is a positive allosteric modulator of GABA<sub>A</sub> receptors.<sup>1</sup> It potentiates GABA-induced currents in LTK cells expressing  $\alpha$ 4 $\beta$ 2 $\gamma$ 2 subunit-containing GABA<sub>A</sub> receptors and CHO cells expressing  $\alpha$ 4 $\beta$ 3 $\delta$  subunit-containing GABA<sub>A</sub> receptors (EC<sub>50</sub>s = 375 and 299 nM, respectively). SAGE-217 (0.3-10 mg/kg, i.p.) reduces the number of seizures induced by pentylenetetrazole (PTZ; Item No. 18682) in mice. It also inhibits electrographic and behavioral seizure activity in a rat model of lithium-pilocarpine induced status epilepticus.

### Reference

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).

#### 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, 10/03/2023

#### 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