

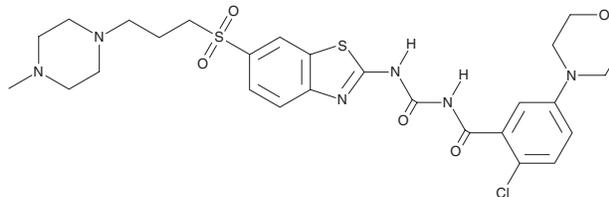
# PRODUCT INFORMATION



## AZ-GHS-22

Item No. 34611

**CAS Registry No.:** 1143020-91-0  
**Formal Name:** 2-chloro-N-[[[6-[[3-(4-methyl-1-piperazinyl)propyl]sulfonyl]-2-benzothiazolyl]amino]carbonyl]-5-(4-morpholinyl)-benzamide  
**MF:** C<sub>27</sub>H<sub>33</sub>ClN<sub>6</sub>O<sub>5</sub>S<sub>2</sub>  
**FW:** 621.2  
**Purity:** ≥98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

AZ-GHS-22 is supplied as a solid. A stock solution may be made by dissolving the AZ-GHS-22 in the solvent of choice. AZ-GHS-22 is soluble in the organic solvent DMSO, which should be purged with an inert gas. The solubility of AZ-GHS-22 in DMSO is approximately 50 mg/ml. AZ-GHS-22 is also soluble in 0.1 M hydrochloride at approximately 25 mg/ml.

### Description

AZ-GHS-22 is an inverse agonist of the growth hormone secretagogue receptor 1a (GHS-R1a), which is also known as the ghrelin receptor.<sup>1</sup> It binds to GHS-R1a with an IC<sub>50</sub> value of 0.77 nM. AZ-GHS-22 decreases food intake in mice by 54% in the first two hours after administration of a 100 mg/kg dose.

### Reference

1. McCoull, W., Barton, P., Brown, A.J.H., *et al.* Identification, optimization, and pharmacology of acylurea GHS-R1a inverse agonists. *J. Med. Chem.* **57**(14), 6128-6140 (2014).

#### 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, 11/22/2022

#### 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