

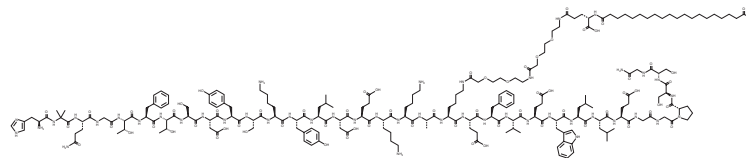
# PRODUCT INFORMATION



## Mazdutide

Item No. 40971

**CAS Registry No.:** 2259884-03-0  
**Formal Name:** L-histidyl-2-methylalanyl-L-glutaminyglycyl-L-threonyl-L-phenylalanyl-L-threonyl-L-seryl-L- $\alpha$ -aspartyl-L-tyrosyl-L-seryl-L-lysyl-L-tyrosyl-L-leucyl-L- $\alpha$ -aspartyl-L- $\alpha$ -glutamyl-L-lysyl-L-lysyl-L-alanyl-N<sup>6</sup>-[N-(19-carboxy-1-oxonadecyl)-L- $\gamma$ -glutamyl-2-[2-(2-aminoethoxy)ethoxy]acetyl-2-[2-(2-aminoethoxy)ethoxy]acetyl]-L-lysyl-L- $\alpha$ -glutamyl-L-phenylalanyl-L-valyl-L- $\alpha$ -glutamyl-L-tryptophyl-L-leucyl-L-leucyl-L- $\alpha$ -glutamylglycylglycyl-L-prolyl-L-seryl-L-seryl-glycinamide  
**Synonyms:** IBI362, LY3305677  
**MF:** C<sub>210</sub>H<sub>322</sub>N<sub>46</sub>O<sub>67</sub>  
**FW:** 4563.1  
**Purity:**  $\geq$ 98%  
**Supplied as:** A 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

Mazdutide is supplied as a solid. Aqueous solutions of mazdutide can be prepared by directly dissolving the solid in aqueous buffers. Mazdutide is sparingly soluble (1-10 mg/ml) in in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

### Description

Mazdutide is a synthetic peptide derivative of the gut peptide hormone oxyntomodulin and a dual agonist of the glucagon-like peptide 1 receptor (GLP-1R) and glucagon receptor (GCGR).<sup>1</sup> It reduces body weight in both *Gcgr*<sup>-/-</sup> and *Glp1r*<sup>-/-</sup> mice.

### Reference

1. Ji, L., Jiang, H., An, P., et al. IBI362 (LY3305677), a weekly-dose GLP-1 and glucagon receptor dual agonist, in Chinese adults with overweight or obesity: A randomised, placebo-controlled, multiple ascending dose phase 1b study. *EClinicalMedicine* **39**, 101088 (2021).

**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, 06/05/2024

### 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