

PRODUCT INFORMATION



ZP3022

Item No. 42636

CAS Registry No.: 1345835-12-2

Formal Name: L-histidylglycyl-L- α -glutamylglycyl-L-threonyl-L-phenylalanyl-L-threonyl-L-seryl-L- α -aspartyl-L-leucyl-L-seryl-L-lysyl-L-glutamyl-L-methionyl-L- α -glutamyl-L- α -glutamyl-L- α -glutamyl-L-alanyl-L-valyl-L-arginyl-L-leucyl-L-phenylalanyl-L-isoleucyl-L- α -glutamyl-L-tryptophyl-L-leucyl-L-lysyl-L-asparaginy-2-[2-(2-aminoethoxy)ethoxy] acetyl-2-[2-(2-aminoethoxy)ethoxy] acetyl-L-tyrosylglycyl-L-tryptophyl-L-leucyl-L- α -aspartyl-L-phenylalaninamide

MF: C₁₉₉H₂₉₄N₄₈O₆₁S

FW: 4,366.9

Purity: $\geq 95\%$

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

ZP3022 is supplied as a solid. A stock solution may be made by dissolving the ZP3022 in the solvent of choice, which should be purged with an inert gas. ZP3022 is soluble (≥ 10 mg/ml) in DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of ZP3022 can be prepared by directly dissolving the solid in aqueous buffers. ZP3022 is slightly soluble (0.1-1 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

ZP3022 is a dual agonist of the glucagon-like peptide 1 receptor (GLP-1R) and cholecystokinin 2 (CCK₂) receptor.¹ It selectively increases cAMP accumulation or ERK phosphorylation in HEK293 cells expressing GLP-1R or the CCK₂ receptor over ERK phosphorylation in HEK293 cells expressing the CCK₁ receptor (EC₅₀s = 0.02, 10.3, and $>1,000$ nM, respectively, for the human receptors). ZP3022 (40 nM) increases the proliferation of rat neonatal β -cells *in vitro*.² It also increases glucose-induced insulin secretion (GSIS) in rat pancreatic islets when used at concentrations of 10 and 100 nM and reduces body weight in rats when administered at a dose of 40 nmol/kg twice per day. ZP3022 (100 nmol/kg per day) increases β -cell mass and total pancreatic mass and improves glucose tolerance in *db/db* mice.¹ It also reduces food intake but not body weight in *db/db* mice.

References

1. Fosgerau, K., Jessen, L., Lind Tolborg, J., *et al.* The novel GLP-1-gastrin dual agonist, ZP3022, increases β -cell mass and prevents diabetes in *db/db* mice. *Diabetes Obes. Metab.* **15**(1), 62-71 (2013).
2. Skarbalienė, J., Rigbolt, K.T., Fosgerau, K., *et al.* In-vitro and in-vivo studies supporting the therapeutic potential of ZP3022 in diabetes. *Eur. J. Pharmacol.* **815**, 181-189 (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, 01/31/2025

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