

PRODUCT INFORMATION



SHR-2042 (sodium salt)

Item No. 42637

Formal Name: N6-(L-histidyl-2-methylalanyl-L- α -glutamylglycyl-L-threonyl-L-phenylalanyl-L-threonyl-L-seryl-L- α -aspartyl-L-valyl-L-seryl-L-seryl-L-tyrosyl-L-leucyl-L- α -glutamyl-L- α -glutamyl-L- α -glutamyl-L-alanyl-L-alanyl-L-lysyl-L- α -glutamyl-L-phenylalanyl-L-isoleucyl-L-alanyl-L-tryptophyl-L-leucyl-L-valyl-L-arginylglycylglycyl-L-prolyl-L-seryl-L-serylglycyl-L-alanyl-L-prolyl-L-prolyl-L-prolyl-L-seryl)-N2-(1-oxohexadecyl)-L-lysine, sodium salt

MF: $C_{205}H_{314}N_{48}O_{63} \cdot XNa$

FW: 4,459.0

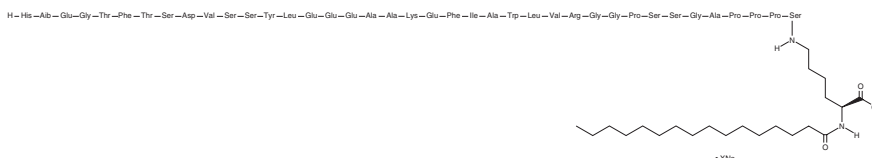
Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years

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



Laboratory Procedures

SHR-2042 (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the SHR-2042 (sodium salt) in the solvent of choice, which should be purged with an inert gas. SHR-2042 (sodium salt) is sparingly soluble (1-10 mg/ml) in DMSO and soluble (≥ 10 mg/ml) in ethanol.

Description

SHR-2042 is an agonist of the glucagon-like peptide 1 receptor (GLP-1R).¹ It is more lipophilic than semaglutide (Item No. 29969). SHR-2042, in combination with sodium N-(8[2-hydroxybenzoyl] amino) caprylate (SNAC; Item No. 36505), has a higher oral bioavailability than semaglutide in monkeys.

Reference

1. Chen, H., Lu, Y., Shi, S., *et al.* Design and development of a new glucagon-like peptide-1 receptor agonist to obtain high oral bioavailability. *Pharm. Res.* **39**(8), 1891-1906 (2022).

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, 04/16/2026

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