

# PRODUCT INFORMATION



## SAR441255 (sodium salt)

Item No. 41175

**Peptide Sequence:** HXHGTFTSDLSKLLK( $\gamma$ Glu-  
 $\gamma$ Glu-palmitic acid)  
EEQRQXEFIEWLKAaGPPSXKPPPK-  
NH<sub>2</sub> (X=2-Aminoisobutyric Acid)

**MF:** C<sub>225</sub>H<sub>355</sub>N<sub>57</sub>O<sub>63</sub> • XNa

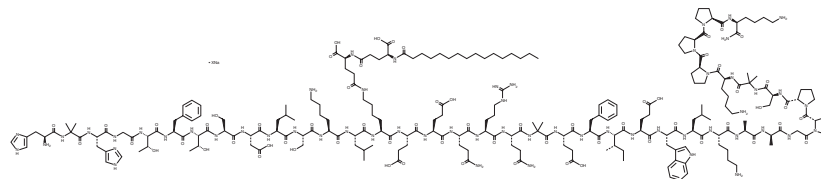
**FW:** 4,866.6

**Purity:**  $\geq$ 95%

**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

SAR441255 (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the SAR441255 (sodium salt) in the solvent of choice, which should be purged with an inert gas. SAR441255 (sodium salt) is slightly soluble (0.1-1 mg/ml) in acetonitrile. It is also slightly soluble (0.1-1 mg/ml) in water.

### Description

SAR441255 is a peptide triple agonist of the glucagon-like peptide 1 receptor (GLP-1R), glucose-dependent insulinotropic polypeptide receptor (GIPR), and glucagon receptor (GCGR).<sup>1</sup> It induces cAMP accumulation in HEK293 cells expressing human GLP-1R, GIPR, or GCGR (EC<sub>50</sub>s = 1.03, 0.73, and 1.01 pM, respectively). SAR441255 (1-30  $\mu$ g/kg) reduces body weight in a mouse model of high-fat diet-induced obesity. It also reduces body weight and blood hemoglobin A1c (HbA1c) levels in high-fat diet-induced obese diabetic cynomolgus monkeys when administered at a dose of 11  $\mu$ g/kg.

### Reference

1. Bossart, M., Wagner, M., Elvert, R., *et al.* Effects on weight loss and glycemic control with SAR441255, a potent unimolecular peptide GLP-1/GIP/GCG receptor triagonist. *Cell Metab.* **34**(1), 59-74 (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, 07/19/2024

#### 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