

# PRODUCT INFORMATION

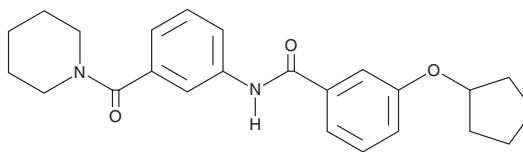


## GLP-1R Modulator L7-028

Item No. 41246

**CAS Registry No.:** 2648317-95-5  
**Formal Name:** N,N,6-triethyl-9,10-didehydro-1-(1-oxopropyl)-ergoline-8 $\beta$ -carboxamide  
**Synonyms:** GLP-1 Receptor Modulator L7-028, Glucagon-like Peptide 1 Receptor Modulator L7-028

**MF:** C<sub>24</sub>H<sub>26</sub>N<sub>2</sub>O<sub>3</sub>  
**FW:** 392.5  
**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

GLP-1R modulator L7-028 is supplied as a solid. A stock solution may be made by dissolving the GLP-1R modulator L7-028 in the solvent of choice, which should be purged with an inert gas. GLP-1R modulator L7-028 is soluble ( $\geq$ 10 mg/ml) in DMSO and sparingly soluble (1-10 mg/ml) in ethanol.

### Description

GLP-1R modulator L7-028 is a positive allosteric modulator of the glucagon-like peptide 1 receptor (GLP-1R).<sup>1</sup> It enhances GLP-1 binding to GLP-1R ( $EC_{50} = 11.01 \mu\text{M}$ ) and cAMP accumulation induced by GLP-1 (7-36) in CHO cells expressing GLP-1R.

### Reference

1. Wang, J., Yang, D., Cheng, X., *et al.* Allosteric modulators enhancing GLP-1 binding to GLP-1R via a transmembrane site. *ACS Chem. Biol.* **16**(11), 2444-2452 (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, 08/27/2024

### 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