

# PRODUCT INFORMATION



## Taspoglutide (acetate)

Item No. 41258

CAS Registry No.: 1022150-16-8

Formal Name: 8-(2-methylalanine)-35-(2-methylalanine)-36-L-argininamide-7-36-glucagon-like peptide I (human), acetate

Synonyms: [Aib<sup>8,35</sup>]hGLP-1(7-36)NH<sub>2</sub>, BIM-51077C

Peptide Sequence: HXEGTFTSDVSSYLEGQAAKEFIAWLVKXR-NH<sub>2</sub> (X = α-aminoisobutyric acid)

MF: C<sub>152</sub>H<sub>232</sub>N<sub>40</sub>O<sub>45</sub> • XC<sub>2</sub>H<sub>4</sub>O<sub>2</sub>

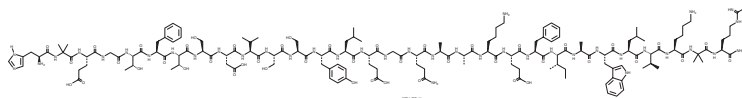
FW: 3,339.7

Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥2 years



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

### Laboratory Procedures

Taspoglutide (acetate) is supplied as a solid. Aqueous solutions of taspoglutide (acetate) can be prepared by directly dissolving the solid in aqueous buffers. Taspoglutide (acetate) 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

Taspoglutide is an agonist of the glucagon-like peptide 1 receptor (GLP-1R; K<sub>i</sub> = 1.1 nM for the human receptor).<sup>1</sup> It induces cAMP accumulation in CHO-K1 cells expressing human GLP-1R (EC<sub>50</sub> = 0.06 nM). Taspoglutide (1 mg/animal per week) decreases blood levels of glucose and increases blood levels of insulin in an oral glucose tolerance test in Zucker diabetic obese rats. It reduces blood levels of gastric inhibitory polypeptide (GIP), plasma levels of triglycerides, and body weight in the same model.<sup>2</sup>

### References

1. Dong, J.Z., Shen, Y., Zhang, J., *et al.* Discovery and characterization of taspoglutide, a novel analogue of human glucagon-like peptide-1, engineered for sustained therapeutic activity in type 2 diabetes. *Diabetes Obes. Metab.* **13**(1), 19-25 (2011).
2. Sebokova, E., Bénardeau, A., Sprecher, U., *et al.* Taspoglutide, a novel human once-weekly analogue of glucagon-like peptide-1, improves glucose homeostasis and body weight in the Zucker diabetic fatty rat. *Diabetes Obes. Metab.* **12**(8), 674-682 (2010).

#### 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, 08/21/2024

#### 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