

PRODUCT INFORMATION



KBP-066 (trifluoroacetate salt)

Item No. 41331

Peptide Sequence: Ac-CSNLSTCXLGRLSQDLH

RLQTYPKTDVGNAP-NH₂

(X=Aminoisobutyric Acid)

MF: C₁₄₇H₂₄₂N₄₆O₄₈S₂ • XCF₃COOH

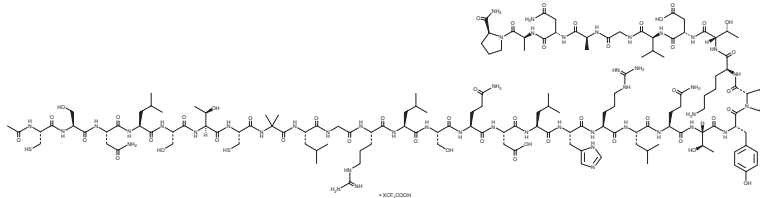
FW: 3,485.9

Purity: ≥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

KBP-066 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the KBP-066 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. KBP-066 (trifluoroacetate salt) 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 KBP-066 (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. KBP-066 (trifluoroacetate salt) 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

KBP-066 is a dual amylin and calcitonin (CT) receptor agonist.¹ It induces β -arrestin recruitment in cells expressing CT receptor subtype a or the amylin 3 receptor (AMY₃) in a concentration-dependent manner. KBP-066 (50 μ g/kg) reduces body weight and food intake in rats fed a high-fat diet. It prevents increases in fasting blood glucose levels in *fa/fa* Zucker diabetic fatty rats.

Reference

1. Sonne, N., Larsen, A.T., Andreassen, K.V., *et al.* The dual amylin and calcitonin receptor agonist, KBP-066, induces an equally potent weight loss across a broad dose range while higher doses may further improve insulin action. *J. Pharmacol. Exp. Ther.* **373**(1), 92-102 (2020).

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, 08/30/2024

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