

PRODUCT INFORMATION



Glucagon Receptor Antagonist 4

Item No. 22164

CAS Registry No.: 1393124-08-7

Formal Name: N-[4-[(1S)-1-[3,5-dimethyl-4-[4-(trifluoromethyl)-1H-pyrazol-1-yl]phenoxy]butyl]benzoyl]-β-alanine

Synonym: GCGR Antagonist IV

MF: $C_{26}H_{28}F_3N_3O_4$

FW: 503.5

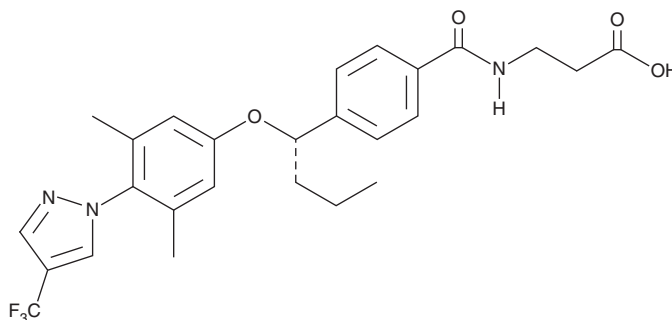
Purity: ≥98%

UV/Vis.: λ_{max} : 234 nm

Supplied as: A crystalline 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

Glucagon receptor antagonist 4 is supplied as a crystalline solid. A stock solution may be made by dissolving the glucagon receptor antagonist 4 in the solvent of choice, which should be purged with an inert gas. Glucagon receptor antagonist 4 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of glucagon receptor antagonist 4 in ethanol is approximately 0.5 mg/ml and approximately 50 mg/ml in DMSO and DMF.

Glucagon receptor antagonist 4 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, glucagon receptor antagonist 4 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Glucagon receptor antagonist 4 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Glucagon receptor antagonist 4 is an antagonist of the glucagon receptor (GCGR; $K_i = 14$ nM for the human recombinant receptor).¹ It is greater than 230-fold selective for GCGR over the glucagon-like peptide 1 (GLP-1) receptor, as well as a panel of 69 receptors, ion channels, uptake sites, and enzymes at 10 μ M. Glucagon receptor antagonist 4 prevents glucagon-induced increases in glucose levels in rat blood when administered at doses of 7.5 and 75 mg/kg.

Reference

1. Guzman-Perez, A., Pfefferkorn, J.A., Lee, E.C., *et al.* The design and synthesis of a potent glucagon receptor antagonist with favorable physicochemical and pharmacokinetic properties as a candidate for the treatment of type 2 diabetes mellitus. *Bioorg. Med. Chem. Lett.* **23**(10), 3051-3058 (2013).

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, 12/16/2022

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