

PRODUCT INFORMATION

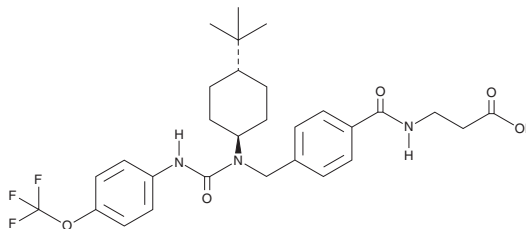


GRA Ex-25

Item No. 29727

CAS Registry No.: 307983-31-9
Formal Name: N-[4-[[[*trans*-4-(1,1-dimethylethyl)cyclohexyl][[4-(trifluoromethoxy)phenyl]amino]carbonyl]amino]methyl]benzoyl]-β-alanine

MF: C₂₉H₃₆F₃N₃O₅
FW: 563.6
Purity: ≥98%
UV/Vis.: λ_{max}: 241 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

GRA Ex-25 is supplied as a crystalline solid. A stock solution may be made by dissolving the GRA Ex-25 in the solvent of choice, which should be purged with an inert gas. GRA Ex-25 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of GRA Ex-25 in ethanol is approximately 10 mg/ml and approximately 25 mg/ml in DMSO and DMF.

GRA Ex-25 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GRA Ex-25 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. GRA Ex-25 has a solubility of approximately 0.1 mg/ml in a 1:8 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

GRA Ex-25 is a glucagon receptor (GCGR) antagonist (IC₅₀s = 55 and 56 nM for the human and rat receptors, respectively).¹ It inhibits adenylate cyclase activity induced by glucagon (Item No. 24204) with a K_i value of 254 nM.² GRA Ex-25 (3 mg/kg, i.v.) reduces glucagon-induced increases in blood glucose levels in rats.¹

References

1. Lau, J., Behrens, C., Sidemann, U.G., *et al.* New β-alanine derivatives are orally available glucagon receptor antagonists. *J. Med. Chem.* **50(1)**, 113-128 (2007)
2. Kurukulasuriya, R., Sorensen, B.K., Link, J.T., *et al.* Biaryl amide glucagon receptor antagonists. *Bioorg. Med. Chem. Lett.* **14(9)**, 2047-2050 (2004).

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, 12/09/2022

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