

PRODUCT INFORMATION



LGD 6972 (ammonium salt)

Item No. 40961

Formal Name: 2-[[4-[(2R)-2-[4'-(1,1-dimethylethyl)[1,1'-biphenyl]-4-yl]-3-oxo-3-[(2',4',6'-trimethyl[1,1'-biphenyl]-4-yl)amino]propyl]benzoyl]amino]-ethanesulfonic acid, monoammonium salt

MF: C₄₃H₄₅N₂O₅S • NH₄

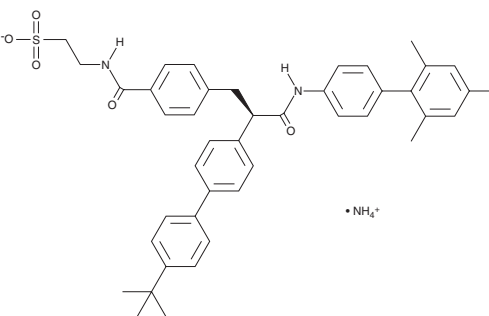
FW: 719.9

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

LGD 6972 (ammonium salt) is supplied as a solid. A stock solution may be made by dissolving the LGD 6972 (ammonium salt) in the solvent of choice, which should be purged with an inert gas. LGD 6972 (ammonium salt) is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

LGD 6972 is a glucagon receptor (GCGR) antagonist.¹ *In vivo*, LGD 6972 reduces glucagon-induced hyperglycemia in wild-type mice and reduces blood glucose levels in mouse models of diabetes.

Reference

1. Vajda, E.G., Logan, D., Lassetter, K., *et al.* Pharmacokinetics and pharmacodynamics of single and multiple doses of the glucagon receptor antagonist LGD-6972 in healthy subjects and subjects with type 2 diabetes mellitus. *Diabetes Obes. Metab.* **19**(1), 24-32 (2017).

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, 11/06/2024

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