

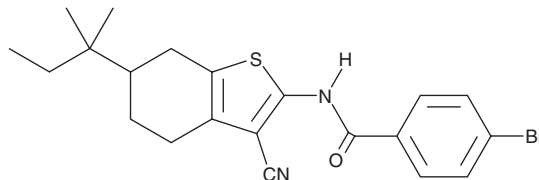
PRODUCT INFORMATION



Glucagon Receptor Antagonist Inactive Control

Item No. 21188

CAS Registry No.: 362482-00-6
Formal Name: 4-bromo-N-[3-cyano-6-(1,1-dimethylpropyl)-4,5,6,7-tetrahydrobenzo[b]thien-2-yl]-benzamide
Synonym: GCGR Antagonist Inactive Control
MF: C₂₁H₂₃BrN₂OS
FW: 431.4
Purity: ≥95%
UV/Vis.: λ_{max}: 246, 322 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 inactive control is supplied as a crystalline solid. A stock solution may be made by dissolving the glucagon receptor antagonist inactive control in the solvent of choice, which should be purged with an inert gas. Glucagon receptor antagonist inactive control is soluble in the organic solvent DMSO.

Description

Glucagon receptor antagonist inactive control is an analog of glucagon receptor antagonist I (Item No. 21223).¹ It inhibits glucagon binding to the glucagon receptor (GCGR) by only 20% in CHO cells expressing human GCGR when used at a concentration of 10 μM. Glucagon receptor antagonist inactive control has no effect on glucagon-induced glycogenolysis in primary human hepatocytes when used at concentrations of 10 or 30 μM.

Reference

1. Qureshi, S.A., Candelore, M.R., Xie, D., *et al.* A novel glucagon receptor antagonist inhibits glucagon-mediated biological effects. *Diabetes* **53**(12), 3267-3273 (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, 06/26/2024

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