

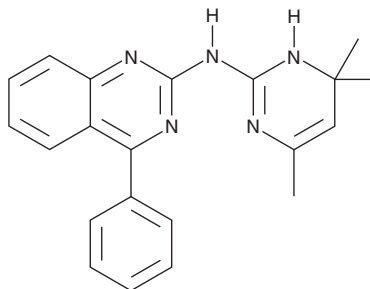
PRODUCT INFORMATION



0990CL

Item No. 45877

CAS Registry No.: 511514-03-7
Formal Name: N-(1,6-dihydro-4,6,6-trimethyl-2-pyrimidinyl)-4-phenyl-2-quinazolinamine
MF: $C_{21}H_{21}N_5$
FW: 343.4
Purity: $\geq 95\%$
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

0990CL is supplied as a solid. A stock solution may be made by dissolving the 0990CL in the solvent of choice, which should be purged with an inert gas. 0990CL is soluble (≥ 10 mg/ml) in ethanol and DMSO.

Description

0990CL is an inhibitor of the G protein subtype $G\alpha_{i1}$.¹ It inhibits $G\alpha_{i1}$ nucleotide exchange by 25% and $G\alpha_q$ nucleotide exchange by $<10\%$ at $300\text{ }\mu\text{M}$. 0990CL ($0.1\text{ }\mu\text{M}$) reverses UK 14304-induced decreases in cAMP in forskolin-stimulated HEK293 cells expressing the α_2 -adrenergic receptor (α_2 -AR).

Reference

1. Appleton, K.M., Bigam, K.J., Lindsey, C.C., *et al.* Development of inhibitors of heterotrimeric $G\alpha_i$ subunits. *Bioorg. Med. Chem.* **22(13)**, 3423-3434 (2015).

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, 06/24/2026

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