

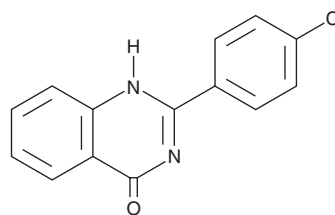
PRODUCT INFORMATION



CQ

Item No. 45825

CAS Registry No.: 7455-77-8
Formal Name: 2-(4-chlorophenyl)-4(3H)-quinazolinone
MF: C₁₄H₉ClN₂O
FW: 256.7
Purity: ≥98%
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

CQ is supplied as a solid. A stock solution may be made by dissolving the CQ in the solvent of choice, which should be purged with an inert gas. CQ is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

CQ is an inhibitor of thermolysin and α -glucosidase (IC_{50} s = 1.25 and 12.5 μ M, respectively).^{1,2}

References

1. Khan, M.T., Khan, R., Wuxiuer, Y., *et al.* Identification of novel quinazolin-4(3H)-ones as inhibitors of thermolysin, the prototype of the M4 family of proteinases. *Bioorg. Med. Chem.* **18(12)**, 4317-4327 (2010).
2. Wei, M., Chai, W.M., Wang, R., *et al.* Quinazolinone derivatives: Synthesis and comparison of inhibitory mechanisms on α -glucosidase. *Bioorg. Med. Chem.* **25(4)**, 1303-1308 (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, 06/29/2026

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