

PRODUCT INFORMATION



C-87

Item No. 36578

CAS Registry No.: 332420-90-3

Formal Name: 4-[2-(4-chloro-3-nitrophenyl)hydrazone]3-phenyl-1-(4-phenyl-2-thiazolyl)-1H-pyrazole-4,5-dione

MF: C₂₄H₁₅ClN₆O₃S

FW: 502.9

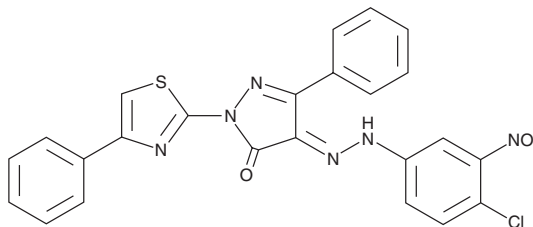
Purity: ≥95%

UV/Vis.: λ_{max}: 247, 403 nm

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

C-87 is supplied as a solid. A stock solution may be made by dissolving the C-87 in the solvent of choice, which should be purged with an inert gas. C-87 is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml (warmed and sonicated).

Description

C-87 is an inhibitor of TNF-α.¹ It binds to TNF-α and reduces the levels of phosphorylated JNK, prevents TNF-α-induced IκBα degradation, and inhibits TNF-α-induced cytotoxicity (IC₅₀ = 8.73 μM) in L-929 cells. It reduces liver inflammation and injury, as well as increases survival, in a mouse model of acute hepatitis induced by LPS and galactosamine (GalN) when administered at a dose of 12.5 mg/kg.

Reference

1. Ma, L., Gong, H., Zhu, H., *et al.* A novel small-molecule tumor necrosis factor α inhibitor attenuates inflammation in a hepatitis mouse model. *J. Biol. Chem.* **289**(18), 12457-12466 (2014).

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, 05/04/2023

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