

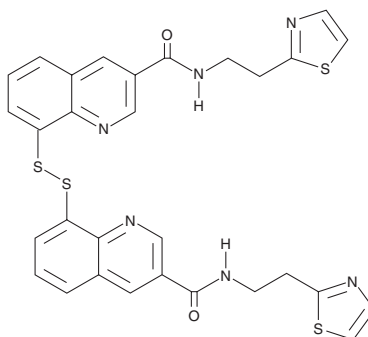
PRODUCT INFORMATION



Capzimin Dimer

Item No. 41364

CAS Registry No.: 2084868-04-0
Formal Name: 8,8'-dithiobis[N-[2-(2-thiazolyl)ethyl]-3-quinolinecarboxamide]
MF: $C_{30}H_{24}N_6O_2S_4$
FW: 628.8
Purity: $\geq 98\%$
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

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

Description

Capzimin dimer is a synthetic intermediate in the synthesis of capzimin, an inhibitor of the proteasome 19S subunit Rpn11.^{1,2}

References

1. Perez, C., Li, J., Parlati, F., *et al.* Discovery of an inhibitor of the proteasome subunit Rpn11. *J. Med. Chem.* **60**(4), 1343-1361 (2017).
2. Li, J., Yakushi, T., Parlati, F., *et al.* Capzimin is a potent and specific inhibitor of proteasome isopeptidase Rpn11. *Nat. Chem. Biol.* **13**(5), 486-493 (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, 08/01/2024

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