

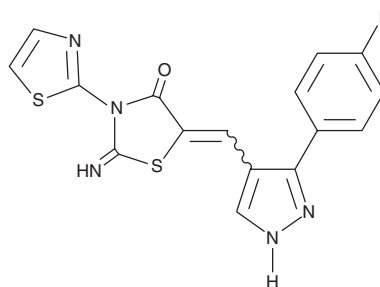
# PRODUCT INFORMATION



## Necrostatin-7

Item No. 10528

**CAS Registry No.:** 351062-08-3  
**Formal Name:** 5-((3-(4-fluorophenyl)-1H-pyrazol-4-yl)methylene)-2-imino-3-(thiazol-2-yl)thiazolidin-4-one  
**Synonym:** Nec-7  
**MF:** C<sub>16</sub>H<sub>10</sub>FN<sub>5</sub>OS<sub>2</sub>  
**FW:** 371.4  
**Purity:** ≥98%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Necrostatin-7 (Nec-7) is supplied as a crystalline solid. A stock solution may be made by dissolving the Nec-7 in the solvent of choice. Nec-7 is soluble in organic solvents such as ethanol and DMSO, which should be purged with an inert gas.

### Description

Necroptosis is a regulated caspase-independent cell death mechanism that results in morphological features resembling necrosis. Serine/threonine kinase activity of the death domain receptor-associated molecule RIP1 is thought to be essential for Fas ligand-induced and tumor necrosis factor- $\alpha$  (TNF- $\alpha$ ) induced necrosis.<sup>1</sup> Nec-7 is a necroptosis inhibitor that is structurally and biologically distinct from necrostatin-1, -3, -4, and -5 as it does not inhibit RIP1 kinase. Nec-7 may target an additional regulatory molecule in this pathway as it inhibits TNF- $\alpha$ -induced necroptosis in a FADD-deficient variant of human Jurkat T cells with an EC<sub>50</sub> value of 10.6  $\mu$ M.<sup>2</sup>

### References

1. Teng, X., Degtarev, A., Jagtap, P., *et al.* Structure-activity relationship study of novel necroptosis inhibitors. *Bioorg. Med. Chem. Lett.* **15**, 5039-5044 (2005).
2. Zheng, W., Degtarev, A., Hsu, E., *et al.* Structure-activity relationship study of a novel necroptosis inhibitor, necrostatin-7. *Bioorg. Med. Chem. Lett.* **18**, 4932-4935 (2008).

#### 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, 10/10/2022

#### 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