

# PRODUCT INFORMATION

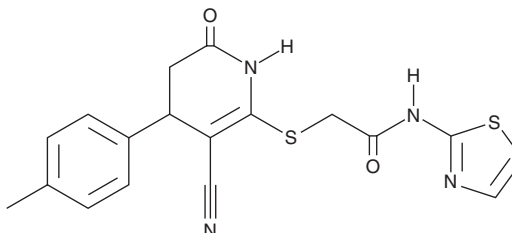


## Necrostatin-34

Item No. 37735

**CAS Registry No.:** 375835-43-1  
**Formal Name:** 2-[[3-cyano-1,4,5,6-tetrahydro-4-(4-methylphenyl)-6-oxo-2-pyridinyl]thio]-N-2-thiazolyl-acetamide

**Synonym:** Nec-34  
**MF:** C<sub>18</sub>H<sub>16</sub>N<sub>4</sub>O<sub>2</sub>S<sub>2</sub>  
**FW:** 384.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 280 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

Necrostatin-34 is supplied as a solid. A stock solution may be made by dissolving the necrostatin-34 in the solvent of choice, which should be purged with an inert gas. Necrostatin-34 is soluble in the organic solvent DMSO.

### Description

Necrostatin-34 is an inhibitor of receptor-interacting serine/threonine kinase 1 (RIPK1; IC<sub>50</sub> = 5.5 μM).<sup>1</sup> It inhibits TNF-α-induced necroptosis in Jurkat cells lacking Fas-associated death domain protein (FADD) and L-929 cells (IC<sub>50</sub>s = 6.67 and 1.35 μM, respectively). Necrostatin-34 (10 μM) inhibits necroptosis induced by the chemical inducer of dimerization for FK506 binding protein (FKBP) fusion proteins AP20187 (Item No. 36808) and the pan-caspase inhibitor Z-VA-DL-D(OMe)-FMK (zVAD-FMK; Item No. 27421) in L-929 cells expressing an FKBP and RIPK1 fusion protein.

### Reference

1. Meng, H., Wu, G., Zhao, X., *et al.* Discovery of a cooperative mode of inhibiting RIPK1 kinase. *Cell Discov.* **7**(1), 41 (2021).

#### 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, 04/19/2023

#### 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