

7-Cl-O-Nec1

Rip1 blocker

Necrostatin-1 analogue with superior potency, selectivity and metabolic stability in blocking RIP1. 7-Cl-O-Nec-1 shows no off-target inhibition of indolamine-2,3-deoxygenase (IDO) in contrast to Necrostatin-1 (Nec-1). 7-Cl-O-Nec-1 showed higher activity in inhibiting necroptosis in Jurkat cells than Necrostatin-1, no non-specific cytotoxicity at high concentrations (100 μ M) and reasonable pharmacokinetic characteristics when used in mice. 7-Cl-O-Nec-1 is recommended for cellular and *in vivo* use over Necrostatin-1.

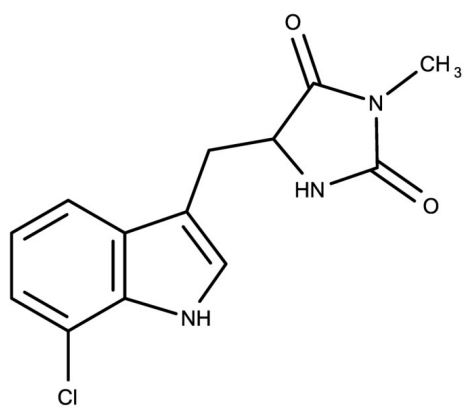
Ordering Information

[Order Online »](#)

| | |
|-----------------|------|
| ENZ-CHM496-0025 | 25mg |
|-----------------|------|

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 5-((7-Chloro-1H-indol-3-yl)methyl)-3-methylimidazolidine-2,4-dione

Appearance Yellow solid.

CAS 852391-15-2

Couple Target RIP

Couple Type Blocker

Formula $C_{13}H_{12}ClN_3O_2$

Identity Determined by NMR.

MW 277.71

Purity ≥98% (TLC)

Solubility Soluble in DMSO (25 mg/ml).

Last modified: February 27, 2025



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)