

# 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  $\mu$ 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.

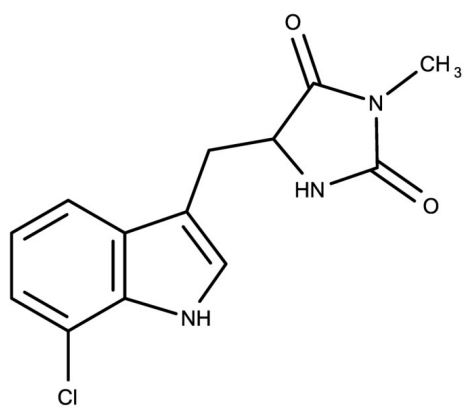
### Ordering Information

[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).



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-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@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)