

7-CL-O-Nec-1

RIPK1 Inhibitor

Catalog # R07-901B

Lot # D3370-1

CAS # 852391-15-2

Product Description

Molecular Formula: C₁₃H₁₂ClN₃O

Appearance: yellow powder

Molecular Weight: 277.71

Purity: >98% (TLC); NMR (Conforms)

Solubilization: May be dissolved DMSO (25 mg/ml)

Alias

Nec-1s; 5-((7-Chloro-1H-indol-3-yl)methyl)-3-methylimidazolidine-2,4-dione

Specific Activity

RIPK1 Inhibitor

Storage and Stability

Store desiccated as supplied at ambient temperature for up to 2 years. Store solutions at -20°C for up to 3 months.

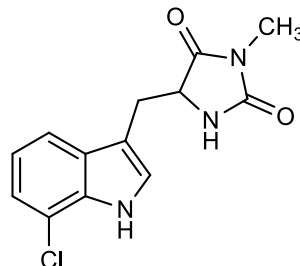
Scientific Background

A necrostatin-1 analog with superior potency², selectivity³ and metabolic stability at RIP1 inhibition (IC₅₀ = 206 vs 494 nM)^{1,2}.

References

1. A Degterev *et al. Nat. Chem. Biol.* 2005 1:112
2. X Teng *et al. Bioorg. Med. Chem. Lett.* 2005 15:5039
3. N Takahashi *et al. Cell Death Dis.* 2012 3:e437

Molecular Structure



7-CL-O-Nec-1

RIPK1 Inhibitor

Catalog Number	R07-901B
CAS Number	852391-15-2
Specific Lot Number	D3370-1
Purity	>98% (TLC); NMR (Conforms)
Format	Yellow Powder
Solubilization	May be dissolved DMSO (25 mg/ml)
Storage & Shipping	Store desiccated as supplied at ambient temperature for up to 2 years. Store solutions at -20°C for up to 3 months

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.