

# 7-CN-7-C-Ino

## Activator of cAMP-independent protein kinase A of Trypanosoma

7-CN-7-C-Ino is a potent and highly specific activator of the cAMP-independent trypanosomal protein kinase A (in *Trypanosoma brucei* EC<sub>50</sub> 6.5 ± 4.6 nM on the purified holoenzyme, KD 8 ± 2 nM to the regulatory subunit PKAR) that has low toxicity *in vivo*. It was also described as a naturally occurring metabolite, called Jaspamycin, isolated from the marine sponge *Jaspis splendens*.

*BLG-C214-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

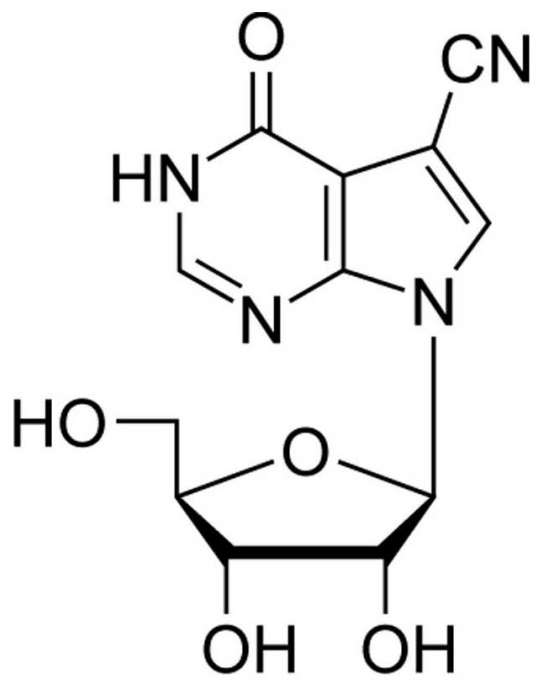
### Ordering Information

[Order Online »](#)

BLG-C214-05	5µmol
BLG-C214-25	5x5µmol

### Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name 7-Cyano-7-deazainosine, Jaspamycin

CAS 22242-96-2

Couple Type Modified nucleotides

Formula  $C_{12}H_{12}N_4O_5$

MW 292.3

Purity > 98% HPLC

Quantity 5 µmol / ~1.5 mg



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-be@enzolifesciences.com](mailto:info-be@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)