

# 8-CEA-Ado

8-CEA-Ado is an analogue of adenosine prepared to be coupled to various structures including proteins. Furthermore, it can be used as a ligand for immobilization to yield affinity gels and is also suitable for conjugation with fluorescent dyes or labels.

*BLG-C092-50 (5 x 10  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

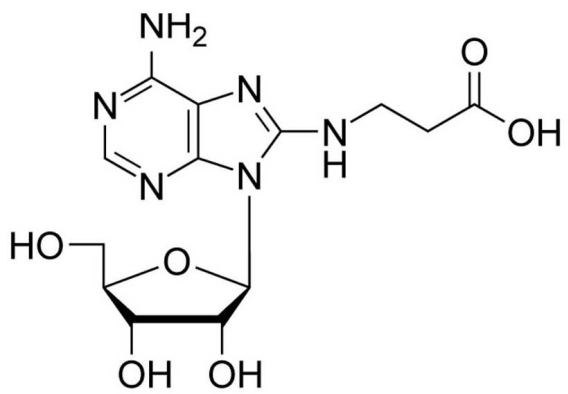
## Ordering Information

[Order Online »](#)

BLG-C092-50	5x10 $\mu$ mol
-------------	----------------

## 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 8-Carboxyethylaminoadenosine . sodium salt

CAS 1151916-53-8

Couple Type Modified nucleotides

Formula  $C_{13}H_{18}N_6O_6$  (free acid)

MW 354.3

Purity > 97% HPLC

Quantity 10µmol≈3.5mg

Solubility Soluble in water (>49mM).

**Technical Info / Product Notes** For the Original Manufacturer's data sheet please [click here](#).



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)