

# 8-Fluo-AET-c-diAMP

## Fluorescent analogue of c-diAMP

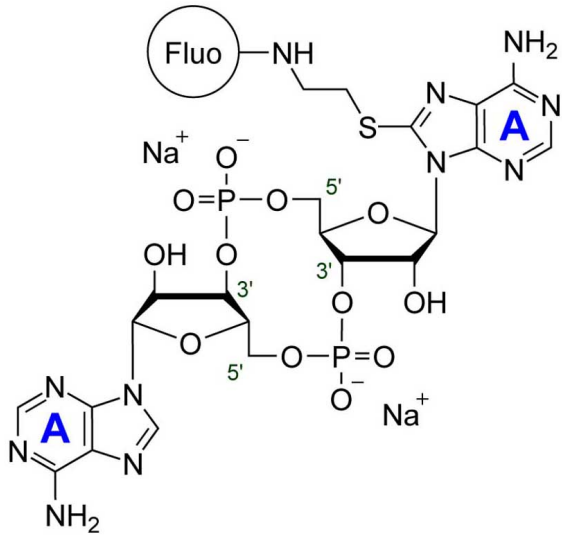
8-Fluo-AET-c-diAMP is a fluorescent analogue of the bacterial second messenger c-diAMP (Prod. No. BLG-C088) with  $\lambda_{exc}$  494 nm and  $\lambda_{em}$  517 nm where one of the two adenine nucleobases is modified by fluoresceine via a short spacer.

*BLG-F014-005 (5 x 0.1  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

### Ordering Information [Order Online »](#)

BLG-F014-001	0.1 $\mu$ mol
BLG-F014-005	5x0.1 $\mu$ mol

### Manuals, SDS & CofA [View Online »](#)



## Handling & Storage

Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 8-(2-[Fluoresceinyl]aminoethylthio)-cyclic diadenosine monophosphate

**Couple Type** Modified nucleotides

**Formula**  $C_{43}H_{39}N_{11}O_{18}P_2S_2$  (free acid)

**MW** 1091.9 (free acid)

**Purity** > 95% HPLC

**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)