

# 2'-Fluo-AHC-c-diGMP

## Fluorescent analogue of c-diGMP

In 2'-Fluo-AHC-c-diGMP, two 5'-GMP units are connected to form a cyclic structure. In addition, 5-carboxyfluorescein has been attached to the 2'-hydroxy group of one ribose via an aminohexylcarbamoyl spacer. 2'-Fluo-AHC-c-diGMP is a fluorescent analogue of the bacterial second messenger c-diGMP (Prod No. [BLG-C057](#)) with  $\lambda_{\text{exc}}$  494 nm and  $\lambda_{\text{em}}$  517 nm.

*BLG-F009-005 (5 x 0.1  $\mu\text{mol}$  pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

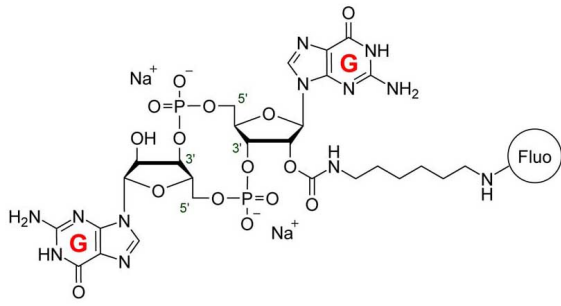
### Ordering Information

[Order Online »](#)

|              |                       |
|--------------|-----------------------|
| BLG-F009-001 | 0.1 $\mu\text{mol}$   |
| BLG-F009-005 | 5x0.1 $\mu\text{mol}$ |

### 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 | 2'-O-(6-[Fluoresceinyl]aminohexylcarbamoyl)-cyclic diguanosine monophosphate . sodium salt |
| CAS              | 1422761-55-4   |
| Couple Type      | Modified nucleotides   |
| Formula          | $C_{48}H_{48}N_{12}O_{21}P_2$ (free acid)  |
| MW               | 1190.9 (free acid)   |
| Purity           | > 95% HPLC   |
| Quantity         | 0.1 $\mu$ mol = ~0.1 mg  |
| Solubility       | Soluble in water (>1.3mM).   |

**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)