

2',2'',2''',2''''',2''''''',2'''''''''-

# Hexa-F-c-hexadAMP

**Analogue of the bacterial second messenger  
c-hexaAMP with increased metabolic  
stability**

2',2'',2''',2''''',2''''''',2'''''''''-Hexa-F-c-hexadAMP is an analogue of the bacterial second messenger c-hexaAMP (Cat. No.: BLG-C332) containing fluoro modifications in all ribose 2'-positions. 2',2'',2''',2''''',2''''''',2'''''''''-Hexa-F-c-hexadAMP was found to be resistant to degradation by the type III CRISPR-associated ribonuclease Csm6 from *Enterococcus italicus* (EiCsm6) while it still activated EiCsm6 (EC50 = 170 nM), albeit more weakly than c-hexaAMP.

*BLG-H038-005 (5 x 0.1 µmol pack size) is not sold in the U.S. or Canada.  
Please [contact us](#) for available options.*

## Ordering Information

[Order Online »](#)

BLG-H038-001	0.1µmol
BLG-H038-005	5x0.1µmol

## 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** 2'-,2"-,2""',2""',2""',2""'-Hexadeoxy-2'-,2""',2""',2""',2""'-hexafluoro-cyclic hexaadenosine monophosphate, sodium salt, cFA6

**Couple Type** Modified nucleotides

**Formula**  $C_{60}H_{66}F_6N_{30}O_{30}P_6$  (free acid)

**MW** 1987.2 (free acid)

**Purity** > 95% HPLC

**Quantity** 0.1 μmol≈~0.2mg



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)