

# 2'-/3'-Fluo-AHC-5'-AMP

## Fluorescent analogue of 5'-AMP

2'-/3'-Fluo-AHC-5'-AMP is an analogue of 5'-AMP in which 5-carboxyfluorescein has been attached either to the ribose 2'-hydroxy group or to the 3'-hydroxy group via an aminohexylcarbamoyl spacer. 2'-/3'-Fluo-AHC-5'-AMP is a fluorescent analogue of 5'-AMP ( $\lambda_{\text{exc}}$  494 nm,  $\lambda_{\text{em}}$  517 nm) and can be used as a calibrator in phosphodiesterase (PDE) assays using 2'-Fluo-AHC-cAMP (FAM-cAMP, Prod. No. [BLG-F003](#)) as a substrate.

*BLG-F010-05 (5 x 1  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

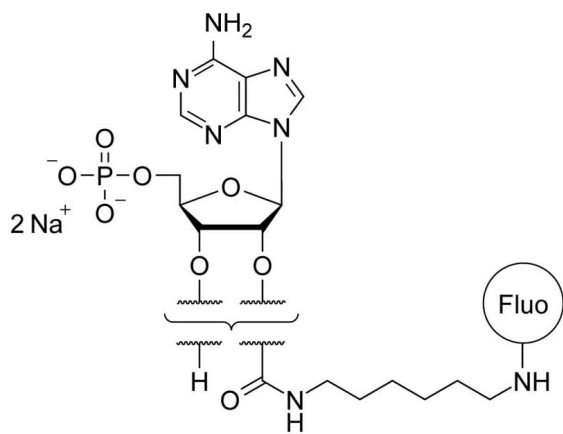
### Ordering Information

[Order Online »](#)

BLG-F010-01	1 $\mu$ mol
BLG-F010-05	5x1 $\mu$ 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	2'-/3'-O-(6-[Fluoresceinyl]aminoethylcarbonyl)adenosine-5'-O-monophosphate . disodium salt, FAM-AMP, FAM-5'-AMP, AMP-Fluorescein conjugate
Couple Type	Modified nucleotides
Formula	$C_{38}H_{38}N_7O_{14}P$ (free acid)
MW	847.7 (free acid)
Purity	> 95% HPLC
Quantity	1 $\mu$ mol = ~0.8 mg
Solubility	Soluble in water (>4mM).
Technical Info / Product Notes	For the Original Manufacturer's data sheet please <a href="#">click here</a> .



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

France  
Phone: +33 472 440 655  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)