

# 8-[DY-547]-AET-cAMP

cyclic AMP conjugated with the fluorescent dye DY-547

Fluorescent cAMP analogue ( $\lambda_{\text{exc}}$  557 nm,  $\lambda_{\text{em}}$  574 nm), e.g. for research focussed on activation studies of HCN2 channels and other cAMP-responsive binding proteins.

*BLG-D109-005 (5 x 0.1  $\mu\text{mol}$  pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

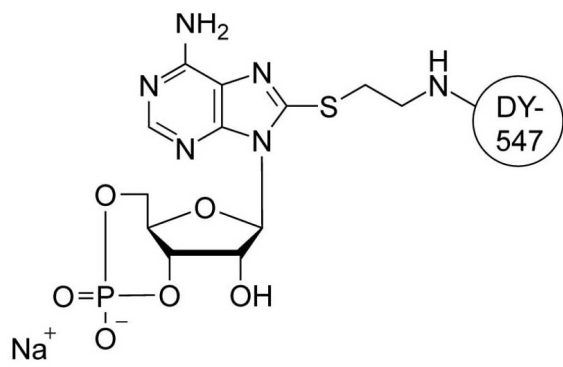
## Ordering Information

[Order Online »](#)

BLG-D109-001	0.1 $\mu\text{mol}$
BLG-D109-005	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	8-(2-[DY-547]-aminoethylthio)adenosine-3',5'-cyclic monophosphate . sodium salt
Couple Type	Modified nucleotides
Formula	$C_{42}H_{49}N_8O_{13}PS_3 \cdot Na_2$
MW	1047.1
Purity	> 98% HPLC
Quantity	0.1µmol≈~0.1mg
Solubility	Soluble in water and aqueous buffers.
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-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)