

# 8-[DY-547P1]-AET-cGMP

cGMP conjugated with the fluorescent dye  
DY-547P1

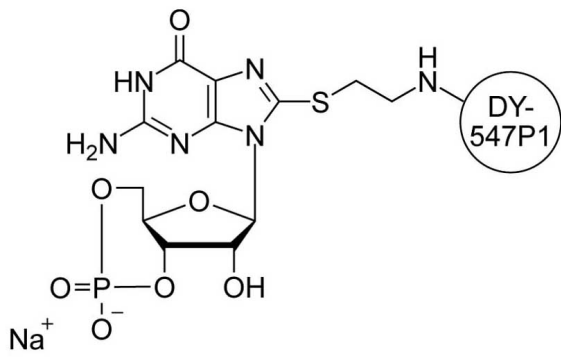
8-[DY-547P1]-AET-cGMP is a fluorescent cGMP analogue ( $\lambda_{exc}$  559 nm,  $\lambda_{em}$  575 nm (EtOH);  $\lambda_{exc}$  551 nm,  $\lambda_{exc}$  565 nm (PBS)) e.g. for research focussed on kinetic studies of CNG ion channels and other cGMP-responsive binding proteins.

*BLG-D224-005 (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-D224-001	0.1 $\mu$ mol
BLG-D224-005	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-[DY-547P1]-aminoethylthio)guanosine-3',5'-cyclic monophosphate . sodium salt
Couple Type	Modified nucleotides
Formula	$C_{44}H_{53}N_8O_{15}PS_3 \cdot Na_2$
MW	1107.1
Purity	> 98% HPLC
Quantity	0.1µmol≈~0.1mg

**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)