

# 2'-TAMRA-AEC-cGMP

cyclic GMP conjugated with the fluorescent dye tetramethylrhodamine

Fluorescent analogue of cGMP with  $\lambda_{\text{exc}}$  555 nm and  $\lambda_{\text{em}}$  580 nm. 2'-TAMRA-AEC-cGMP can be used as a substrate for phosphodiesterases (PDEs) in corresponding assays. This type of fluorescent analogue is also available with a longer 9-atom spacer (2'-TAMRA-AHC-cGMP, Prod. No. BLG-T017).

*BLG-T018-025 (5 x 0.5  $\mu\text{mol}$  pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

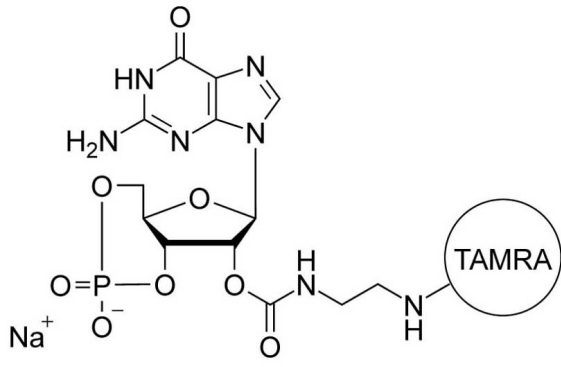
## Ordering Information

[Order Online »](#)

BLG-T018-005	0.5 $\mu\text{mol}$
BLG-T018-025	5x0.5 $\mu\text{mol}$

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Handling	Protect from light.
Long Term Storage	-80°C
Shipping	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	2'-O-(2-[Tetramethylrhodaminy]aminoethylcarbamoyl)guanosine-3',5'-cyclic monophosphate . sodium salt
Couple Type	Modified nucleotides
Formula	$C_{38}H_{37}N_9O_{12}P \cdot Na$
MW	865.7
Purity	> 95% HPLC
Quantity	0.5µmol≈433 µg
Solubility	Soluble in water (≥5mM).
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)