

2'-TAMRA-AHC-cGMP

cyclic GMP conjugated with the fluorescent dye tetramethylrhodamine

Fluorescent analogue of cGMP with λ_{exc} 555 nm and λ_{em} 580 nm. 2'-TAMRA-AHC-cGMP can be used as a substrate for phosphodiesterases (PDEs) in corresponding assays. This type of fluorescent analogue is also available with a shorter 5-atom spacer (2'-TAMRA-AEC-cGMP, Prod. No. BLG-T018).

BLG-T017-025 (5 x 0.5 μmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

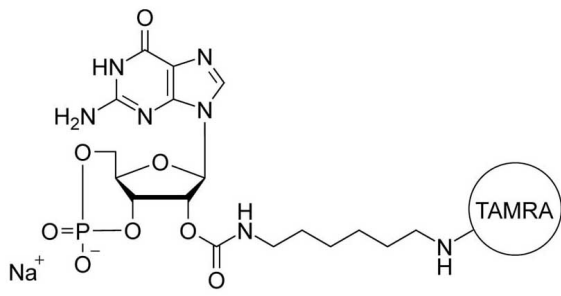
Ordering Information

[Order Online »](#)

BLG-T017-005	0.5 μmol
BLG-T017-025	5x0.5 μmol

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-(6-[Tetramethylrhodaminy]aminohexylcarbamoyl)guanosine-3',5'-cyclic monophosphate . sodium salt
Couple Type	Modified nucleotides
Formula	$C_{42}H_{45}N_9O_{12}P \cdot Na$
MW	921.8
Purity	> 95% HPLC
Quantity	0.5µmol≈461 µg
Solubility	Soluble in water (≥2mM).
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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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