

2'-/3'-TAMRA-AEC-5'- GMP

Fluorescent analogue of 5'-GMP

Fluorescent analogue of guanosine-5'-monophosphate (λ_{exc} 555 nm; λ_{em} 580 nm).

*BLG-T020-025 (5 x 0.5 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

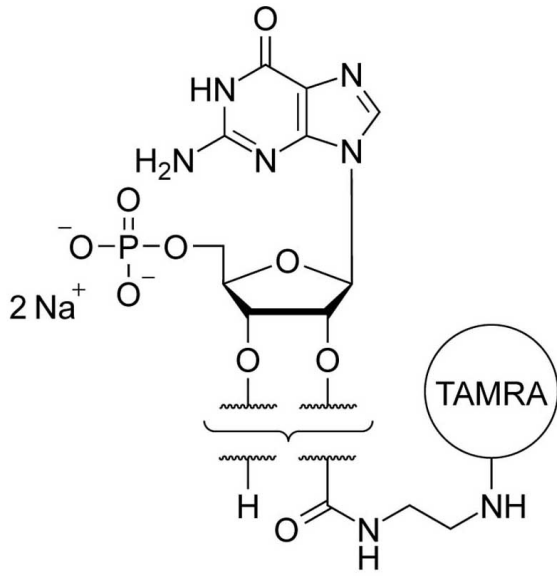
Ordering Information

[Order Online »](#)

BLG-T020-005	0.5 μ mol
BLG-T020-025	5x0.5 μ 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'-/3'-O-(6-[Tetramethylrhodaminyl]aminoethylcarbamoyl)guanosine-5'-O-monophosphate . disodium salt
Formula	$C_{38}H_{40}N_9O_{13}P$ (free acid)
MW	861.8
Purity	> 95% HPLC



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