

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

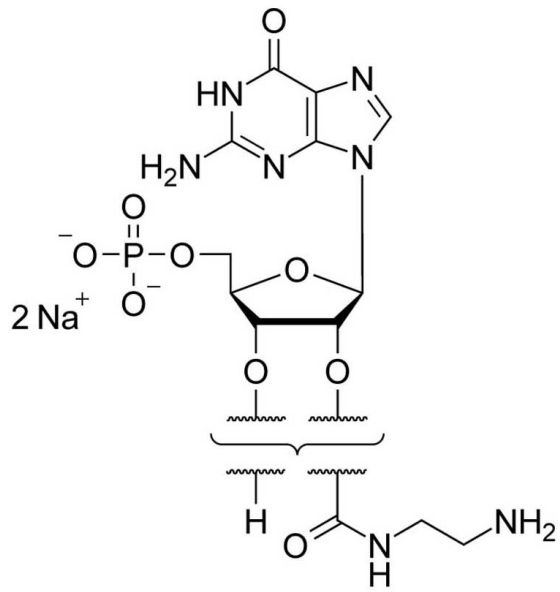
Analogue of guanosine-5'-O-monophosphate, useful as ligand for immobilization and for coupling of various dyes and labels. Detailed technical information available. This ligand is also available with a longer spacer (2'-/3'-AHC-5'-GMP; Cat. No. A 174).

*BLG-A170-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

Ordering Information [Order Online »](#)

BLG-A170-05	5µmol
BLG-A170-25	5x5µmol

Manuals, SDS & CofA [View Online »](#)



Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 2'-/3'-O-(2-Aminoethylcarbamoyl)guanosine-5'-O-monophosphate . disodium salt

Couple Type Modified nucleotides

Formula $C_{13}H_{20}N_7O_9P$

MW 449.3

Purity > 97% 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