

2-NH₂-6-Cl-5'-PuMP

Functionalized 5'-GMP for modification of position 6

BLG-A084-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.

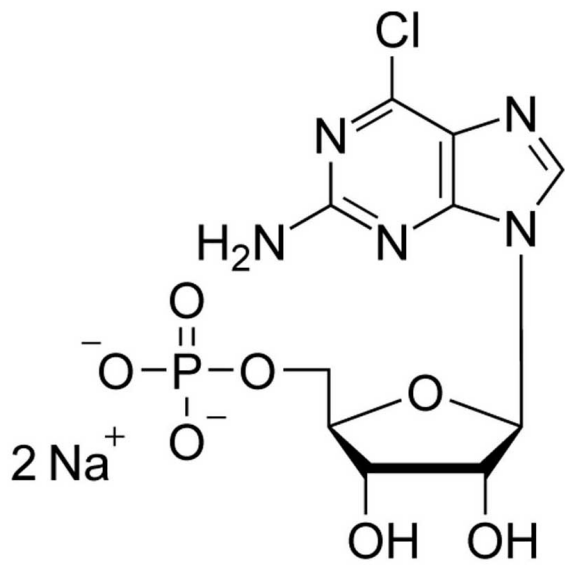
Ordering Information

[Order Online »](#)

| | |
|-------------|---------|
| BLG-A084-05 | 5µmol |
| BLG-A084-25 | 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-Amino-6-chloropurine riboside-5'-O-monophosphate . disodium salt

CAS 16321-98-5

Couple Type Modified nucleotides

Formula $C_{10}H_{13}ClN_5O_7P$

MW 381.7

Purity > 97% HPLC

Quantity 5µmol≈1.9mg

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