

6-Cl-5'-PuMP. Na

Reactive precursor of 6- modified 5'-AMP analogues

Analogue of 5'-AMP with reactive chlorine for introduction of modifications in position 6 of the adenine nucleobase.

BLG-C032-50 (5 x 10 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

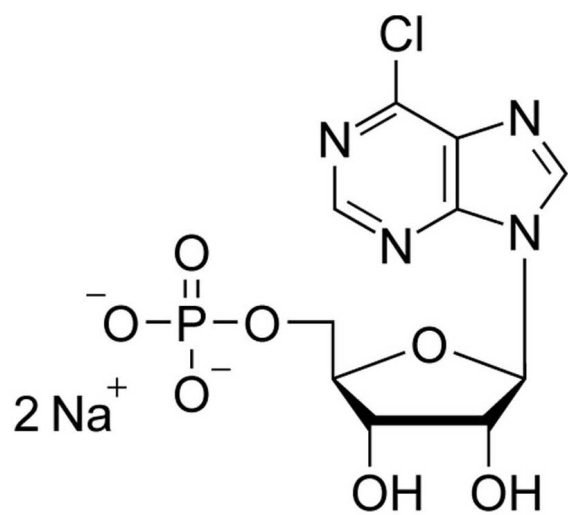
Ordering Information

[Order Online »](#)

BLG-C032-10	10µmol
BLG-C032-50	5x10µmol

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 6-Chloropurine riboside-5'-O-monophosphate . disodium salt

CAS 5843-59-4

Couple Type Modified nucleotides

Formula $C_{10}H_{12}ClN_4O_7P$ (free acid)

MW 366.7 (free acid)

Purity > 97% HPLC

Quantity 10µmol≈3.7mg

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