

6-Bnz-5'-AMP

Lipophilic analogue of adenosine-5'-O-monophosphate and potential metabolite of N6-Benzoyl-cAMP (Cat. No. B 009). For the corresponding nucleoside or nucleobase please inquire. 6-Bnz-5'-AMP shows some cytokinin activity in tobacco and soybean tissue culture assays

*BLG-B109-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

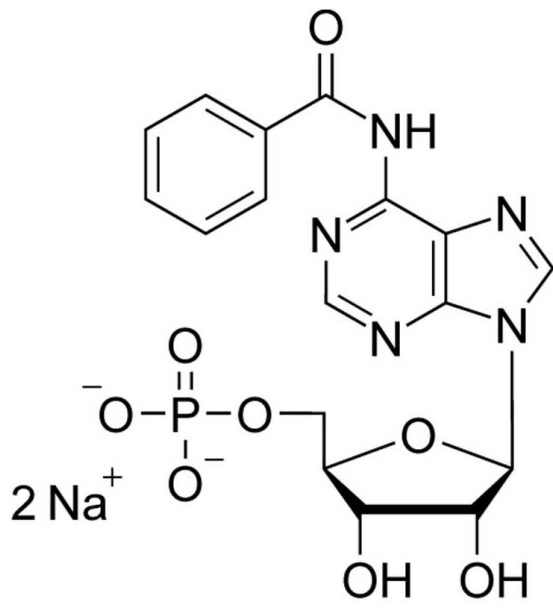
Ordering Information

[Order Online »](#)

BLG-B109-05	5 μ mol
BLG-B109-25	5x5 μ mol

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 N6-Benzoyladenine-5'-O-monophosphate . disodium salt

CAS 40871-55-4

Couple Type Modified nucleotides

Formula $C_{17}H_{18}N_5O_8P$ (free acid)

MW 451.3

Purity > 97% HPLC

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