

7-CH-5'-AMP

Potential substrate, competitive inhibitor or regulator of enzymes that interact with adenosine- 5'- monophosphate.

BLG-D032-50 (5 x 10 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

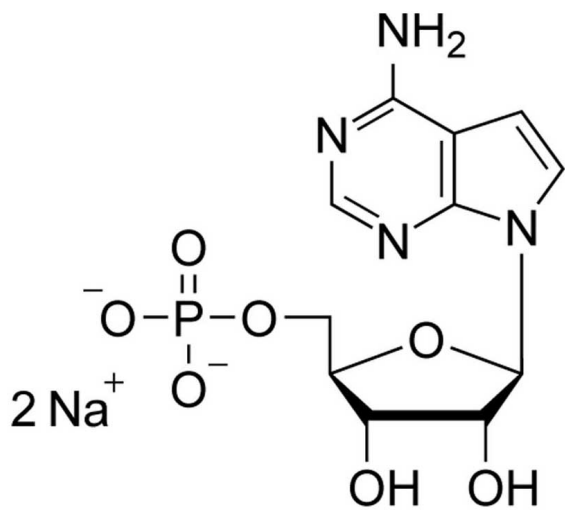
Ordering Information

[Order Online »](#)

BLG-D032-50	5x10 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 7-Deazaadenosine-5'-O-monophosphate . disodium salt, 5'-TuMP, Tubercidin-5'-O-monophosphate

CAS 16719-46-3

Couple Type Modified nucleotides

Formula $C_{11}H_{15}N_4O_7P$ (free acid)

MW 346.2 (free acid)

Purity > 97% HPLC

Quantity 10µmol≈3.5mg

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