

2-MeS-Ado

2-MeS-Ado is an inhibitor of human betaine-homocysteine methyltransferase expression and is also suitable as starting material for 2-methylthio nucleotide analogues. The corresponding 5'-monophosphate 2-MeS-5'-AMP (Prod. No. [BLG-M031](#)) and the 5'-diphosphate 2-MeS-ADP (Prod. No. [BLG-M020](#)) as well as the 5'-triphosphate 2-MeS-ATP (Prod. No. [BLG-M021](#)) are available as well.

BLG-M033-50 (5 x 10 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

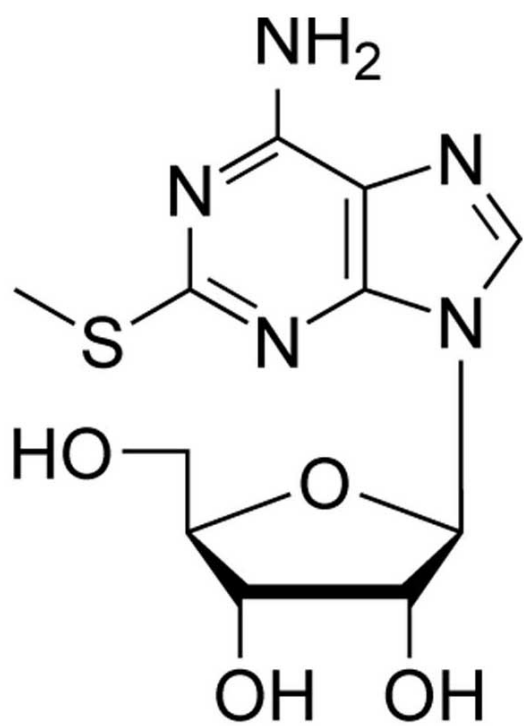
Ordering Information

[Order Online »](#)

BLG-M033-50	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 2-Methylthioadenosine

CAS 4105-39-9

Couple Type Modified nucleotides

Formula $C_{11}H_{15}N_5O_4S$

MW 313.4

Purity > 98% HPLC

Quantity 10µmol≈3.1mg

Solubility Soluble in DMSO or DMF.

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