

2'-O-MS-cAMP

Cyclic AMP analogue, suitable for coupling of various ligands including proteins for ELISA.

BLG-M011-125 (5 x 25 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

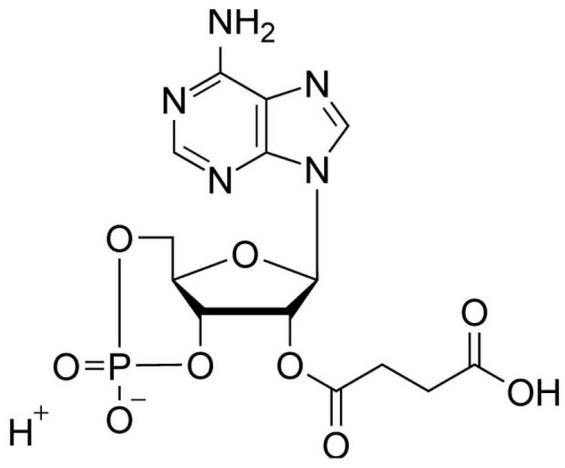
Ordering Information

[Order Online »](#)

BLG-M011-25	25µmol
BLG-M011-125	5x25µmol

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'-Monosuccinyladenosine-3',5'-cyclic monophosphate

CAS 36940-87-1

Couple Type Modified nucleotides

Formula $C_{14}H_{16}N_5O_9P$

MW 429.3

Purity > 98% HPLC

Quantity 25µmol≈11mg

Solubility The free acid form of 2'-O-MS-cAMP is soluble in organic solvents such as DMSO (16.7mM) and DMF (3.3mM), or in aqueous buffer at pH 6-7 (100mM Na₂HPO₄, pH 7.0, 25mM), leading to the respective salt form of 2'-O-MS-cAMP. Concentrations have been tested at ambient temperature and can be considered as minimum concentrations usually obtainable.

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