

m^{2,2,7}G/TMG

m^{2,2,7}G is an analogue of guanosine in which the amino hydrogens in position 2 and in position 7 of the nucleobase are replaced by methyl groups. m^{2,2,7}G can be found in the cap structures at the 5' ends of snRNAs and, far less often, of mRNAs. It can be used as a reference for analysis of methylated snRNAs or mRNAs.

BLG-T009-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

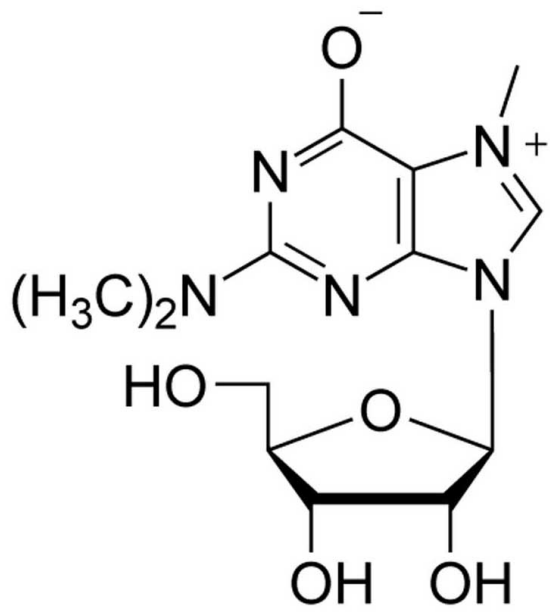
Ordering Information

[Order Online »](#)

BLG-T009-05	5µmol
BLG-T009-25	5x5µmol

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 N2,N2,7-Trimethylguanosine (internal salt)

CAS 40027-70-1

Couple Type Modified nucleotides

Formula $C_{13}H_{19}N_5O_5$ (internal salt)

MW 325.3 (internal salt)

Purity > 95% HPLC

Quantity 5µmol≈1.6mg

Reconstitution Soluble in water or buffers. Please rinse tube walls carefully and preferably use ultrasonic or vortex to achieve total and uniform mixing.

Solubility Soluble in water or buffers.

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