

BCMCM-caged 8-Br-cGMP

Equatorial isomer. BCMCM-caged 8-Br-cGMP is an only weakly fluorescent, caged precursor of 8-bromo cyclic cGMP, which upon irradiation, releases both 8-Br-cGMP and a fluorescent coumarin analogue.

BLG-B019-005 (5 x 0.1 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

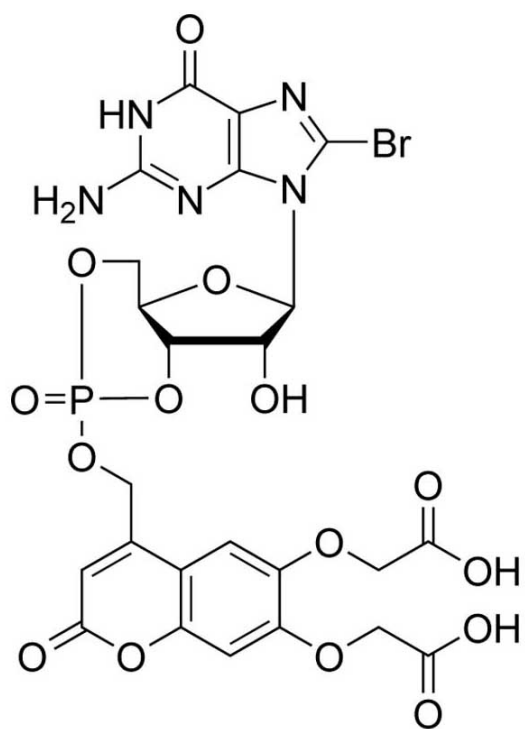
Ordering Information

[Order Online »](#)

BLG-B019-001	0.1 μ mol
BLG-B019-005	5x0.1 μ mol

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Handling	Protect from light.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	[6,7-Bis(carboxymethoxy)coumarin-4-yl]methyl-8-bromoguanosine-3',5'-cyclic monophosphate
Couple Type	Modified nucleotides
Formula	$C_{24}H_{21}BrN_5O_{15}P \cdot 3H_2O$
MW	730.4 . 54.0
Purity	>98% by HPLC
Quantity	0.1µmol≈~78mg
Solubility	Excellent solubility 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](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)