

DEACM-caged cGMP

Caged cyclic GMP, releasing a fluorescent coumarin upon uncaging

Axial isomer. DEACM-caged cGMP is an only weakly fluorescent, caged precursor of the second messenger cyclic GMP, which upon irradiation, releases both cGMP and a fluorescent coumarin analogue. High lipophilicity and membrane permeability; but still reasonable solubility in aqueous systems.

BLG-D043-005 (5 x 0.1 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

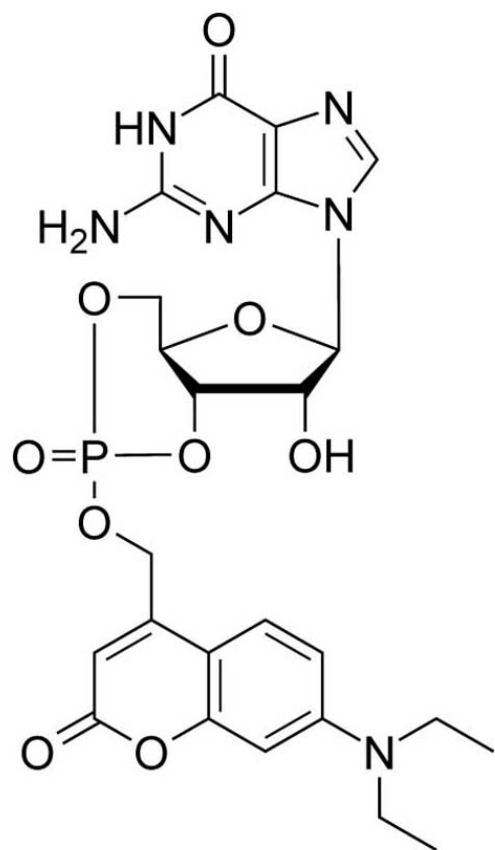
Ordering Information

[Order Online »](#)

BLG-D043-001	0.1 μ mol
BLG-D043-005	5x0.1 μ mol

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name (7-Diethylaminocoumarin-4-yl)methyl guanosine-3',5'-cyclic monophosphate

CAS 339291-42-8

Couple Type Modified nucleotides

Formula $C_{24}H_{27}N_6O_9P$

MW 628.5

Purity > 98% HPLC

Quantity 0.1 μmol ≈ 63 μg

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)