

DMACM-caged 8-Br-cGMP

Caged 8-Br-cGMP releasing fluorescent coumarin upon irradiation

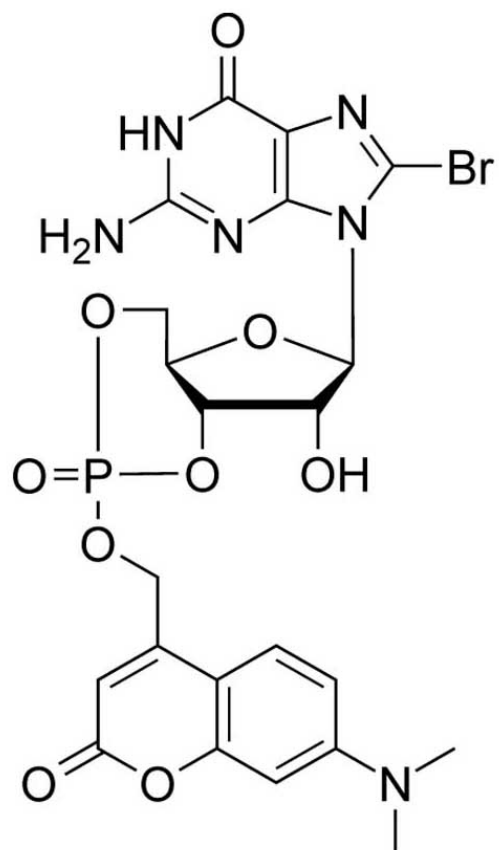
Axial isomer. DMACM-caged cAMP is an only weakly fluorescent, caged precursor of 8- bromo cyclic GMP, which upon irradiation, releases both 8-Br-cGMP and a fluorescent coumarin analogue. High lipophilicity and membrane permeability; thus only limited solubility in aqueous systems.

BLG-D045-005 (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-D045-001	0.1 μ mol
BLG-D045-005	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-Dimethylaminocoumarin-4-yl)methyl-8-bromoguanosine-3',5'-cyclic monophosphate

Formula $C_{22}H_{22}BrN_6O_9P$

MW 634.3

Purity > 98% HPLC

Quantity 0.1 μmol ≈ 60 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

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