

# 8-Br-2',3'-cGMP

Membrane-permeable analogue of 2',3'-cGMP

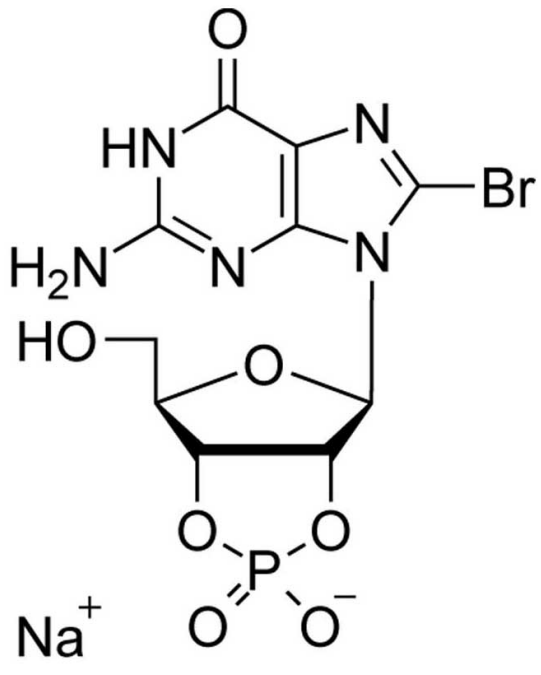
8-Br-2',3'-cGMP is considered to be a membrane-permeable analogue of 2',3'-cGMP (Biolog Cat. No. G 025). 2',3'-cGMP is a positional isomer of the natural second messenger 3',5'-cGMP (Biolog Cat. No. G 001) and was found to exist in mammals under physiological conditions

*BLG-B292-25 (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-B292-05	5µmol
BLG-B292-25	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** 8-Bromoguanosine-2',3'-cyclic monophosphate . sodium salt

**CAS** 52216-45-2

**Couple Type** Modified nucleotides

**Formula**  $C_{10}H_{10}BrN_5O_7P \cdot Na$

**MW** 446.09

**Purity** > 97% HPLC

**Quantity** 5µmol are equivalent to ~2.2 mg

**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)