

8-Fluo-AET-c-diGMP

Fluorescent analogue of c-diGMP

8-Fluo-AET-c-diGMP is a fluorescent analogue of the bacterial second messenger c-diGMP (Cat. No. C 057) with λ_{exc} 494 nm and λ_{em} 517 nm where one of the two guanine nucleobases is modified by fluoresceine via a short spacer.

BLG-F019-005 (5 x 0.1 μmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

Ordering Information

[Order Online »](#)

BLG-F019-001	0.1 μmol
BLG-F019-005	5x0.1 μmol

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-(2-[Fluoresceinyl]aminoethylthio)-cyclic diguanosine monophosphate . sodium salt

Appearance Crystallized or lyophilized solid

Couple Type Modified nucleotides

Formula $C_{43}H_{39}N_{11}O_{20}P_2S$ (free acid)

MW 1123.9

Purity > 95% HPLC

Quantity 0.1µmol are equivalent to ~0.1 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

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