

8-Fluo-cAMP

cyclic AMP conjugated with fluorescein

Fluorescent cAMP analogue. Excitation 494nm, Emission 517nm.
Potent activator of cAMP-dependent protein kinases and expected to be for cAMP-gated ion channels as well.

*BLG-F002-05 (5 x 1 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

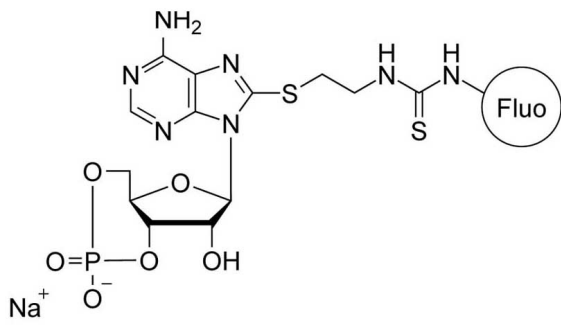
Ordering Information

[Order Online »](#)

BLG-F002-01	1 μ mol
BLG-F002-05	5x1 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	8-[[2-[(Fluoresceinylthioureido)amino]ethyl]thio]adenosine-3',5'-cyclic monophosphate . sodium salt
Appearance	Crystallized or lyophilized sodium salt.
CAS	293296-57-8
Couple Type	Modified nucleotides
Formula	$C_{33}H_{27}N_7O_{11}PS_2 \cdot Na$
MW	815.7
Purity	> 95% HPLC
Quantity	1µmol≈0.8mg
Solubility	Soluble in water, methanol, PBS (pH 7.4) and other 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

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