

8-[AF488]-AHA-cAMP

Fluorescent analogue of cAMP

Fluorescent cAMP analogue (λ_{exc} 493 nm, λ_{em} 517 nm).

*BLG-A412-005 (5 x 0.1 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

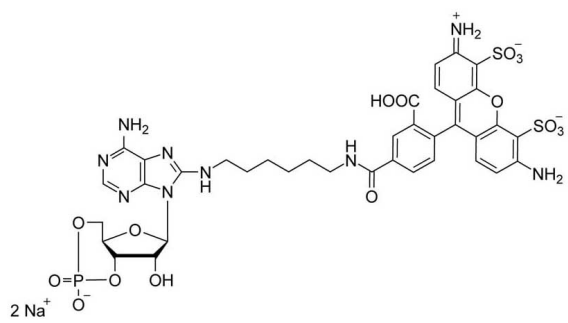
Ordering Information

[Order Online »](#)

BLG-A412-001	0.1 μ mol
BLG-A412-005	5x0.1 μ mol

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-(6-[AF488]-aminohexylamino)adenosine-3',5'-cyclic monophosphate, sodium salt, Alexa Fluor® 488 cAMP, F-cAMP

Couple Type Modified nucleotides

Formula $C_{37}H_{36}N_9O_{16}PS_2 \cdot Na_2$

MW 1003.8

Purity > 98% HPLC

Quantity 0.1 µmol ≈ 0.1 mg



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