

8-[(6-Amino)hexyl]- amino-cAMP – Cy5

Ordering Information

[Order Online »](#)

JBS-NU-851-CY5	160µl
----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 8-[(6-Amino)hexyl]-amino-adenosine-3',5'-cyclic monophosphate, labeled with Cy5, Triethylammonium salt

Appearance blue solution

Couple Target cAMP

Couple Type Modified nucleotides

Emission Maximum 670 nm

Excitation Maximum 649 nm

Formula $C_{49}H_{64}N_9O_{13}PS_2$ (free acid)

Formulation solution in water

MW 1082.19 (free acid)

Purity ≥ 95 % (HPLC)

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-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-de@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-uk@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com