

8-[DY-547]-AET-2'-O-Me-cAMP

Fluorescent Epac agonist

*BLG-D089-005 (5 x 0.1 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

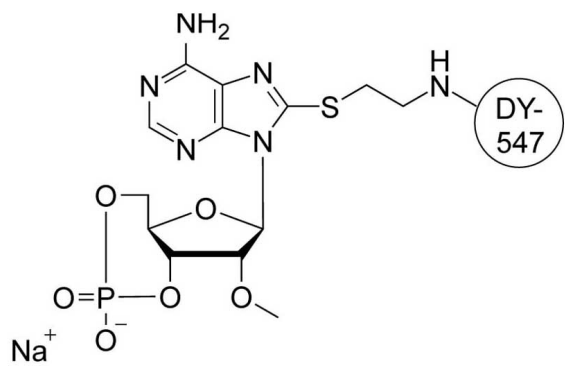
Ordering Information

[Order Online »](#)

| | |
|--------------|-----------------|
| BLG-D089-001 | 0.1 μ mol |
| BLG-D089-005 | 5x0.1 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-[DY-547]-aminoethylthio)-2'-O-methyladenosine-3',5'-cyclic monophosphate . sodium salt |
| Couple Type | Modified nucleotides |
| Formula | $C_{43}H_{51}N_8O_{13}PS_3 \cdot Na_2$ |
| MW | 1061.1 |
| Purity | > 98% HPLC |
| Quantity | 0.1 μmol ≈ 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