

Br-ANT-dATP

Fluorescent dATP analogue

Analogue of 2'-deoxyadenosine-5'-triphosphate with only moderate intrinsic fluorescence (λ_{exc} 345 nm, λ_{em} ~425 nm). The fluorophore is supposed to have a certain sensitivity for its environment and can change its spectral properties upon binding. Useful for research into dATP-dependent receptor proteins. Potential inhibitor of various adenylyl cyclase isoforms and soluble guanylyl cyclase (sGC).

BLG-B145-05 (5 x 1 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

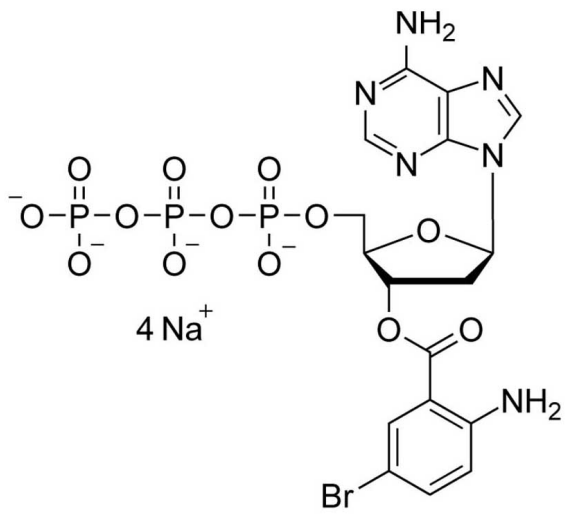
Ordering Information

[Order Online »](#)

BLG-B145-01	1 μ mol
BLG-B145-05	5x1 μ mol

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 3'-O-(5-Bromoanthraniloyl)-2'-deoxyadenosine-5'-O-triphosphate . sodium salt

Couple Type Modified nucleotides

Formula $C_{17}H_{20}BrN_6O_{13}P_3$ (free acid)

Formulation 10 mM aqueous solution

MW 689.2 (free acid)

Purity > 95% HPLC

Quantity 1 μ mol is equivalent to ~0.7 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