

2',2'',2'''-Tri-F-c-tridAMP

Fluorinated version of the bacterial second messenger c-triAMP

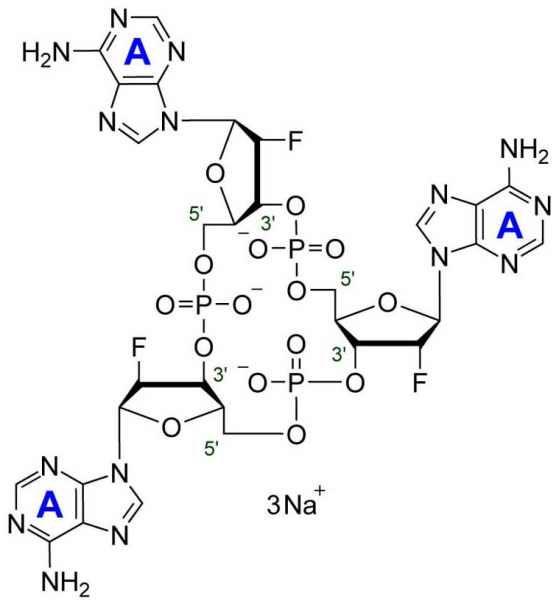
2',2'',2'''-Tri-F-c-tridAMP is an analogue of the bacterial second messenger c-triAMP (Cat. No.: BLG-C362) containing fluoro modifications in all ribose 2'-positions which are expected to make the compound resistant to degradation by nucleases.

BLG-T064-005 (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-T064-001	0.1 μ mol
BLG-T064-005	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 2'-,2"-,2"-Trideoxy-2'-,2",2"-trifluoro-cyclic triadenosine monophosphate . sodium salt, Trifluoro-c-triAMP

Appearance Crystallized or lyophilized solid.

Couple Type Modified nucleotides

Excitation Maximum 259 nm

Formula $C_{30}H_{33}F_3N_{15}O_{15}P_3$ (free acid)

MW 993.6 (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-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