

Ap₄C

Ap₄C is a dinucleoside polyphosphate containing both, a purine nucleobase as well as a pyrimidine nucleobase moiety. Ap₄C exhibits agonist activity at the human P2Y1 and P2Y2 receptors (Shaver et al. 2005).

BLG-A178-025 (5 x 0.5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

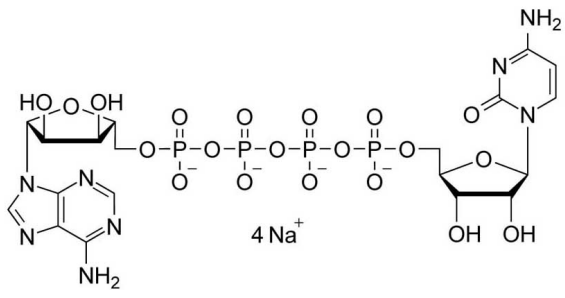
Ordering Information

[Order Online »](#)

BLG-A178-005	0.5µmol
BLG-A178-025	5x0.5µmol

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 P1-(5'-Adenosyl)-P4-(5'-cytidyl)-tetraphosphate . tetrasodium salt, Cp4A

CAS 1352921-13-1

Formula $C_{19}H_{28}N_8O_{20}P_4$ (free acid)

MW 812.4

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