

PRODUCT INFORMATION



Cyclic ADP-Ribose (ammonium salt)

Item No. 21417

Formal Name: 1-β-D-ribofuranosyl-adenosine
5'-(trihydrogen diphosphate),
intramol. P',5''-ester, diammonium salt

Synonyms: cADP-Ribose, cADPR

MF: C₁₅H₁₉N₅O₁₃P₂ • 2NH₄

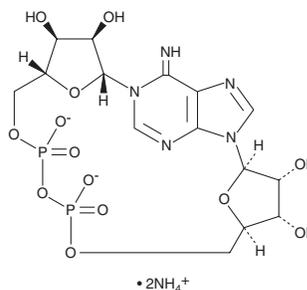
FW: 575.4

Purity: ≥98%

Supplied as: A solid

Storage: -80°C

Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Cyclic ADP-Ribose (cADP-ribose) (ammonium salt) is supplied as a solid. A stock solution may be made by dissolving the cADP-ribose (ammonium salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

cADP-Ribose is an endogenous metabolite of NAD⁺ that mobilizes the release of stored Ca²⁺ in the endoplasmic reticulum *via* ryanodine receptors in various cell types.¹⁻⁵ This second messenger is generated *via* the cADP-ribose synthases CD38 and CD157.^{3,5,6} cADP-Ribose may also trigger the cell surface Ca²⁺ influx channel TRPM2 in a temperature-dependent manner.⁷ *In vitro*, cADP-ribose modulates Ca²⁺ signaling in rat and mouse cardiomyocytes treated with isoproterenol (Item No. 15592), and treatment with this metabolite at 100 μM under heat stress conditions induces the release of oxytocin (Item No. 11799) from the mouse hypothalamus.^{4,6}

References

1. Lee, H.C., Walseth, T.F., Bratt, G.T., *et al.* Structural determination of a cyclic metabolite of NAD⁺ with intracellular Ca²⁺-mobilizing activity. *J. Biol. Chem.* **264**(3), 1608-1615 (1989).
2. Guse, A.H. Biochemistry, biology, and pharmacology of cyclic adenosine diphosphoribose (cADPR). *Curr. Med. Chem.* **11**, 847-855 (2004).
3. Houtkooper, R.H., Cantó, C., Wanders, R.J., *et al.* The secret life of NAD⁺: An old metabolite controlling new metabolic signaling pathways. *Endocr. Rev.* **31**(2), 194-223 (2010).
4. Zhong, J., Amina, S., Liang, M., *et al.* Cyclic ADP-ribose and heat regulate oxytocin release *via* CD38 and TRPM2 in the hypothalamus during social or psychological stress in mice. *Front. Neurosci.* **10**(304), 1-14 (2016).
5. Nikiforov, A., Kulikova, V., and Ziegler, M. The human NAD metabolome: Functions, metabolism and compartmentalization. *Crit. Rev. Biochem. Mol. Biol.* **50**(4), 284-297 (2015).
6. Gul, R., Park, D.-R., Shawl, A.I., *et al.* Nicotinic acid adenine dinucleotide phosphate (NAADP) and cyclic ADP-ribose (cADPR) mediate Ca²⁺ signaling in cardiac hypertrophy induced by β-adrenergic stimulation. *PLoS One* **11**(3), e0149125 (2016).
7. Lee, H.C. Cyclic ADP-ribose and nicotinic acid adenine dinucleotide phosphate (NAADP) as messengers for calcium mobilization. *J. Biol. Chem.* **287**(38), 31633-31640 (2012).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 04/05/2018

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM