

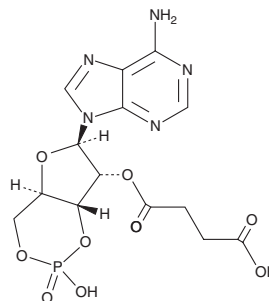
PRODUCT INFORMATION



2'-O-Monosuccinyladenosine-3',5'-cyclic monophosphate

Item No. 30535

CAS Registry No.:	36940-87-1
Formal Name:	cyclic 3',5'-(hydrogen phosphate) 2'-(hydrogen butanedioate) adenosine
Synonyms:	2'-O-Succinyl-cAMP, 2'-O-Succinyl-3',5'-cyclic AMP
MF:	C ₁₄ H ₁₆ N ₅ O ₉ P
FW:	429.3
Purity:	≥98%
UV/Vis.:	λ _{max} : 258 nm
Supplied as:	A crystalline solid
Storage:	-20°C
Stability:	≥2 years



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

Laboratory Procedures

2'-O-Monosuccinyladenosine-3',5'-cyclic monophosphate is supplied as a crystalline solid. Aqueous solutions of 2'-O-monosuccinyladenosine-3',5'-cyclic monophosphate can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 2'-O-monosuccinyladenosine-3',5'-cyclic monophosphate in 1M NH₄OH is approximately 50 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

2'-O-Monosuccinyladenosine-3',5'-cyclic monophosphate is an immunogenic derivative of cAMP that has been used to generate anti-cAMP antisera and antibodies.^{1,2}

References

1. Ezan, E., Mamas, S., Rougeot, C., *et al.* Strategies for developing specific and sensitive hapten radioimmunoassays. *Radioimmunoassay in Basic and Clinical Pharmacology*. Patrono, C. and Peskar, B.A., editors, 1st edition, Springer (1987).
2. Denizot, F.C., Hirn, M.H., and Delaage, M.A. Relationship between affinity and kinetics in a hapten-antibody reaction. Studies on anti-cyclic AMP antibodies. *Mol. Immunol.* **16(7)**, 509-513 (1979).

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, 05/20/2020

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