

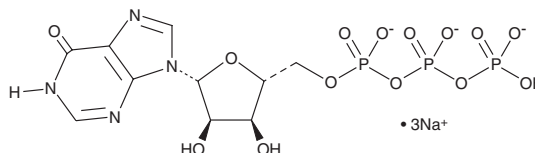
PRODUCT INFORMATION



Inosine 5'-triphosphate (sodium salt)

Item No. 23063

CAS Registry No.: 35908-31-7
Formal Name: inosine 5'-(tetrahydrogen triphosphate), trisodium salt
Synonyms: ITP, NSC 20266
MF: $C_{10}H_{12}N_4O_{14}P_3 \cdot 3Na$
FW: 574.1
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 249 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

Inosine 5'-triphosphate (ITP) (sodium salt) is supplied as a crystalline solid. Aqueous solutions of ITP (sodium salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of ITP (sodium salt) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

ITP is a nucleoside triphosphate that functions as an alternative substrate for ATPases and GTPases.^{1,2} It has been used to study activation and binding kinetics of nucleoside interaction with various ATPases and GTPases.

References

1. Hanson, R.L. and Kennedy, E.P. Energy-transducing adenosine triphosphatase from *Escherichia coli*: Purification, properties, and inhibition by antibody. *J. Bacteriol.* **114**(2), 772-781 (1973).
2. Pieper, U., Schweitzer, T., Groll, D.H., et al. The GTP-binding domain of McrB: More than just a variation on a common theme? *J. Mol. Biol.* **292**(3), 547-556 (1999).

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, 12/05/2022

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