

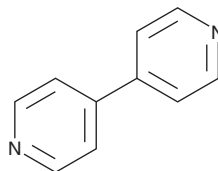
PRODUCT INFORMATION



4,4'-Bipyridyl

Item No. 45661

CAS Registry No.: 553-26-4
Synonyms: BPY, 4,4'B, 4,4'-Dipyridyl, NSC 404423
MF: $C_{10}H_8N_2$
FW: 156.2
Purity: $\geq 98\%$
Supplied as: A 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

4,4'-Bipyridyl is supplied as a solid. A stock solution may be made by dissolving the 4,4'-bipyridyl in the solvent of choice, which should be purged with an inert gas. 4,4'-Bipyridyl is soluble (≥ 10 mg/ml) in ethanol and DMSO.

Description

4,4'-Bipyridyl is a building block that has been used as a ligand in coordination polymer chemistry and the synthesis of metal-organic frameworks (MOFs).^{1,2} It is also a synthetic intermediate that has been used in the synthesis of the herbicide paraquat (Item No. 38521).³

References

1. Lu, W.-G., Jiang, L., Feng, X.-L., *et al.* Three 3D coordination polymers constructed by Cd(II) and Zn (II) with imidazole-4,5-dicarboxylate and 4,4'-bipyridyl building blocks. *Cryst. Growth Des.* **6(2)**, (2006).
2. Yamazui, D., Uchida, K., Koyama, S., *et al.* Syntheses, structures, and properties of coordination polymers with 2,5-dihydroxy-1,4-benzoquinone and 4,4'-bipyridyl synthesized by in situ hydrolysis method. *ACS Omega* **7(22)**, 18259-18266 (2022).
3. Michel-Buono, M., Buono, J.P., Serre, G., *et al.* In vitro cytotoxic effects of 4,4'-bipyridyl on normal human keratinocytes. *Cell Biol. Toxicol.* **13(3)**, 193-204 (1997).

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, 05/12/2026

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