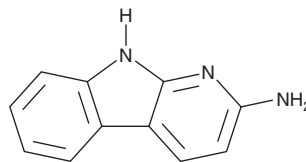


PRODUCT INFORMATION

2-Amino-9H-pyrido[2,3-b]indole

Item No. 42114

CAS Registry No.: 26148-68-5
Formal Name: 9H-pyrido[2,3-b]indol-2-amine
Synonyms: AaC, 2-Amino- α -Carboline
MF: C₁₁H₉N₃
FW: 183.2
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

2-Amino-9H-pyrido[2,3-b]indole (AaC) is supplied as a solid. A stock solution may be made by dissolving the AaC in the solvent of choice, which should be purged with an inert gas. AaC is soluble (≥ 10 mg/ml) in DMSO.

Description

AaC is a mutagenic heterocyclic amine that has been found in cooked foods.¹ It is mutagenic to *S. typhimurium* in the Ames test and induces sister chromatid exchanges (SCEs) in NL3 lymphoblastoid cells.^{1,2} Dietary administration of AaC (0.08%) induces hepatic and blood vessel tumor formation in mice.³ It has also been found in the urine of cigarette smokers.⁴

References

1. Yoshida, D., Matsumoto, T., and Nishigata, H. Effect of heating methods on mutagenic activity and yield of mutagenic compounds in pyrolysis products of protein. *Arg. Biol. Chem.* **44**(2), 253-255 (1979).
2. Tohda, H., Oikawa, A., Kawachi, T., et al. Induction of sister-chromatid exchanges by mutagens from amino acid and protein pyrolysates. *Mutat. Res.* **77**(1), 65-69 (1980).
3. Ohgaki, H., Matsukura, N., Morino, K., et al. Carcinogenicity in mice of mutagenic compounds from glutamic acid and soybean globulin pyrolysates. *Carcinogenesis* **5**(6), 815-819 (1984).
4. Turesky, R.J., Yuan, J.M., Wang, R., et al. Tobacco smoking and urinary levels of 2-amino-9H-pyrido[2,3-b]indole in men of Shanghai, China. *Cancer Epidemiol. Biomarkers Prev.* **16**(8), 1554-1560 (2007).

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, 11/04/2024

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