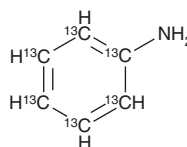


# PRODUCT INFORMATION



Aniline- $^{13}\text{C}_6$   
Item No. 30523

CAS Registry No.: 100849-37-4  
Formal Name: benzenamine-1,2,3,4,5,6- $^{13}\text{C}_6$   
Synonyms: Aminobenzene- $^{13}\text{C}_6$ , Benzenamine- $^{13}\text{C}_6$ ,  
Phenylamine- $^{13}\text{C}_6$   
MF:  $[\text{C}_6\text{H}_7\text{N}]$   
FW: 99.1  
Purity:  $\geq 98\%$   
UV/Vis.:  $\lambda_{\text{max}}$ : 290 nm  
Supplied as: A neat oil  
Storage:  $-20^\circ\text{C}$   
Stability:  $\geq 4$  years



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

## Laboratory Procedures

Aniline- $^{13}\text{C}_6$  is supplied as a neat oil. A stock solution may be made by dissolving the aniline- $^{13}\text{C}_6$  in the solvent of choice, which should be purged with an inert gas. Aniline- $^{13}\text{C}_6$  is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of aniline- $^{13}\text{C}_6$  in these solvents is approximately 30 mg/ml.

Aniline- $^{13}\text{C}_6$  is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, aniline- $^{13}\text{C}_6$  should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Aniline- $^{13}\text{C}_6$  has a solubility of approximately 0.2 mg/ml in a 1:4 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

Aniline- $^{13}\text{C}_6$  is a building block.<sup>1,2</sup> It has been used in the synthesis of  $^{13}\text{C}_6$ -labeled derivatives of tryptophan and indole, as well as  $^{13}\text{C}_6$ -labeled benzimidazoles.

## References

1. Ilić, N. and Cohen, J.D. Synthesis of  $[\text{C}_6\text{H}_7\text{N}]$ -isotopomers of indole and tryptophan for use in the analysis of indole-3-acetic acid biosynthesis. *J. Label. Compd. Radiopharm.* **47(10)**, 635-646 (2004).
2. Donahue, M.G., Jentsch, N.G., and Simons, C.R. Synthesis of  $[\text{C}_6\text{H}_7\text{N}]$ 3,4-diaminobenzoic acid as a precursor for stable isotope labeled benzimidazoles. *Tetrahedron Lett.* **58(17)**, 1692-1694 (2017).

### 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, 09/28/2022

## 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