PRODUCT INFORMATION



Sulfanilic Acid

Item No. 33451

CAS Registry No.:	121-57-3	
Formal Name:	4-amino-benzenesulfonic acid	
Synonyms:	4-Aminobenzenesulfonic Acid, p-Aminobenzenesulfonic Acid, NSC 7170, 4-Sulfoaniline, p-Sulfoaniline	O \\OH S\\
MF:	C ₆ H ₇ NO ₃ S	
FW:	173.2	
Purity:	≥95%	H ₂ N
Supplied as:	A crystalline 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

Sulfanilic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the sulfanilic acid in the solvent of choice, which should be purged with an inert gas. Sulfanilic acid is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of sulfanilic acid in these solvents is approximately 1 mg/ml.

Sulfanilic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, sulfanilic acid should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Sulfanilic acid has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Sulfanilic acid is a synthetic intermediate.^{1,2} It has been used in the synthesis of activated carbon for biodiesel preparation, as well as various nanocomposites.

References

- 1. Malins, K., Kampars, V., Brinks, J., et al. Synthesis of activated carbon based heterogenous acid catalyst for biodiesel preparation. Appl. Catal. B-Environ. 176-177, 553-558 (2015).
- 2. Zhang, L., Shi, T., Wu, S., et al. Sulfonated graphene oxide: The new and effective material for synthesis of polystyrene-based nanocomposites. Colloid Polym. Sci. 291(9), 2061-2068 (2013).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFFTY 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, 12/06/2022

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