

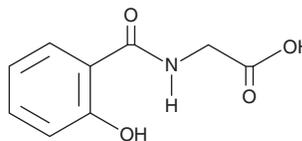
PRODUCT INFORMATION



Salicyluric Acid

Item No. 34813

CAS Registry No.: 487-54-7
Formal Name: N-(2-hydroxybenzoyl)-glycine
Synonyms: 2-Hydroxyhippuric Acid, NSC 524135, N-Salicyloylglycine
MF: C₉H₉NO₄
FW: 195.2
Purity: ≥95%
UV/Vis.: λ_{max}: 237 nm
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

Salicyluric acid is supplied as a solid. A stock solution may be made by dissolving the salicyluric acid in the solvent of choice, which should be purged with an inert gas. Salicyluric acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of salicyluric acid in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Salicyluric acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, salicyluric acid should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Salicyluric acid has a solubility of approximately 0.09 mg/ml in a 1:10 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Salicyluric acid is a major metabolite of both aspirin (Item No. 70260) and salicylic acid.^{1,2}

References

1. Finnie, M.D., Ersser, R.S., Seakins, J.W., *et al.* The occurrence and identification of o-hydroxyhippuric acid (salicyluric acid) in the urine of sick children. *Clin. Chim. Acta* **70(1)**, 171-178 (1976).
2. Needs, C.J. and Brooks, P.M. Clinical pharmacokinetics of the salicylates. *Clin. Pharmacokinet.* **10(2)**, 164-177 (1985).

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, 10/11/2022

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