

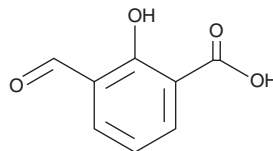
PRODUCT INFORMATION



3-Formylsalicylic Acid

Item No. 23368

CAS Registry No.: 610-04-8
Formal Name: 3-formyl-2-hydroxy-benzoic acid
Synonym: NSC 55753
MF: C₈H₆O₄
FW: 166.1
Purity: ≥95%
UV/Vis.: λ_{max}: 221, 329 nm
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.

Description

3-Formylsalicylic acid is a non-competitive inhibitor of estrone sulfatase (IC₅₀ = 150 nM; K_i = 120 nM), the enzyme that converts estrone sulfate to estrone.¹ 3-Formylsalicylic acid has been used in the synthesis of Schiff bases.^{2,3}

References

1. Jütten, P., Schumann, W., Härtl, A., *et al.* Thiosemicarbazones of formyl benzoic acids as novel potent inhibitors of estrone sulfatase. *J. Med. Chem.* **50(15)**, 3661-3666 (2007).
2. Sarkar, S. and Dey, K. Synthesis and spectroscopic characterization of some transition metal complexes of a new hexadentate N₂S₂O₂ Schiff base ligand. *Spectrochim. Acta. A. Mol. Biomol. Spectrosc.* **62(1-3)**, 383-393 (2005).
3. Hadi, J.S. and Althahabi, N.K. Synthesis, spectroscopic characterization and biological activity of some new sulfa drug Schiff bases. *Res. J. Pharm. Biol. Chem. Sci.* **5(3)**, 856-866 (2014).

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, 11/17/2022

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