

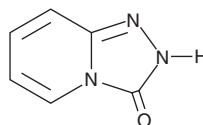
PRODUCT INFORMATION



Triazolopyridinone

Item No. 34393

CAS Registry No.: 6969-71-7
Formal Name: 1,2,4-triazolo[4,3-a]pyridin-3(2H)-one
Synonym: NSC 68462
MF: $C_6H_5N_3O$
FW: 135.1
Purity: $\geq 98\%$
UV/Vis.: λ_{\max} : 217 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

Triazolopyridinone is supplied as a solid. A stock solution may be made by dissolving the triazolopyridinone in the solvent of choice, which should be purged with an inert gas. Triazolopyridinone is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of triazolopyridinone in these solvents is approximately 25 and 15 mg/ml, respectively.

Triazolopyridinone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, triazolopyridinone should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Triazolopyridinone has a solubility of approximately 0.14 mg/ml in a 1:6 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Triazolopyridinone is an intermediate in the synthesis of the antidepressant trazadone.¹ It is also a potential impurity found in commercial preparations of trazadone.²

References

1. Fan, G., Liu, H., Fan, B., *et al.* Trazodone as an efficient corrosion inhibitor for carbon steel in acidic and neutral chloride-containing media: Facile synthesis, experimental and theoretical evaluations. *J. Mol. Liquids* **311**, 113302 (2020).
2. Pai, N.R. and Pusalkar, D.A. Development and validation of liquid chromatographic method for Trazodone hydrochloride. *J. Chem. Pharm. Res.* **2(2)**, 478-488 (2010).

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, 11/22/2022

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