

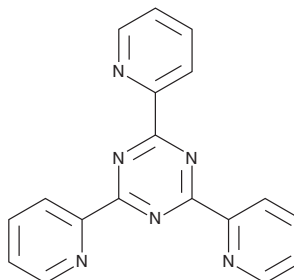
PRODUCT INFORMATION



TPTZ

Item No. 28274

CAS Registry No.: 3682-35-7
Formal Name: 2,4,6-tri-2-pyridinyl-1,3,5-triazine
Synonyms: NSC 112125, 2,4,6-Tripyridyl-s-triazine
MF: C₁₈H₁₂N₆
FW: 312.3
Purity: ≥95%
UV/Vis.: λ_{max}: 245, 284 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.

Laboratory Procedures

TPTZ is supplied as a crystalline solid. A stock solution may be made by dissolving the TPTZ in the solvent of choice, which should be purged with an inert gas. TPTZ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of TPTZ in ethanol is approximately 20 mg/ml and approximately 10 mg/ml in DMSO and DMF.

TPTZ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TPTZ should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. TPTZ 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

TPTZ is a colorimetric reagent commonly used in the detection of iron.^{1,2} TPTZ forms a complex with Fe(II) that can be quantified by colorimetric detection at 594 nm as a measure of iron concentration.

References

1. Collins, P.F. and Diehl, H. 2,4,6-Tripyridyl-s-triazine as a reagent for iron. *Anal. Chem.* **31(11)**, 1862-1867 (1959).
2. Dehghani, Z., Dadfarnia, S., Shabani, A.M.H., *et al.* Determination of iron species by combination of solvent assisted-dispersive solid phase extraction and spectrophotometry. *Anal. Bioanal. Chem. Res.* **2(1)**, 13-21 (2015).

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, 12/14/2022

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