

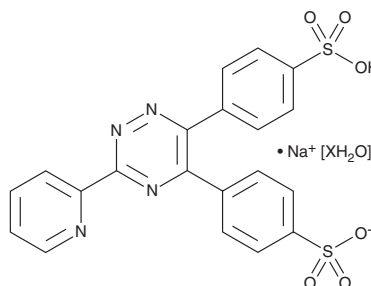
PRODUCT INFORMATION



Ferrozine (hydrate)

Item No. 31465

CAS Registry No.: 1266615-85-3
Formal Name: 4,4'-[3-(2-pyridinyl)-1,2,4-triazine-5,6-diyl]bis-benzenesulfonic acid, monosodium salt, hydrate
MF: C₂₀H₁₃N₄O₆S₂ • Na [XH₂O]
FW: 492.5
Purity: ≥98%
UV/Vis.: λ_{max}: 287 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

Ferrozine (hydrate) is supplied as a crystalline solid. Aqueous solutions of ferrozine (hydrate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of ferrozine (hydrate) in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Ferrozine is a colorimetric reagent commonly used in the detection of iron.^{1,2} Ferrozine forms a complex with ferrous iron that can be quantified by colorimetric detection at 562 nm as a measure of iron concentration.

References

1. Stookey, L.L. Ferrozine-A new spectrophotometric reagent for iron. *Anal. Chem.* **42(7)**, 779-781 (1970).
2. Jeitner, T.M. Optimized ferrozine-based assay for dissolved iron. *Anal. Biochem.* **454**, 36-37 (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, 11/28/2022

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