

PRODUCT INFORMATION

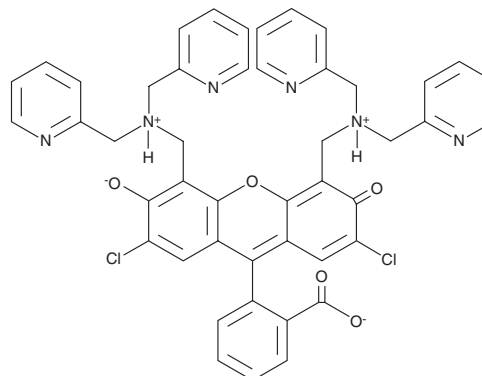


Zinpyr-1

Item No. 15122

CAS Registry No.: 288574-78-7
Formal Name: 4',5'-bis[[bis(2-pyridinylmethyl)amino]methyl]-2',7'-dichloro-3',6'-dihydroxy-spiro[isobenzofuran-1(3H),9'-[9H]xanthen]-3-one

Synonyms: ZP-1
MF: C₄₆H₃₆Cl₂N₆O₅
FW: 823.7
Purity: ≥90%
UV/Vis.: λ_{max}: 232, 513 nm
Ex./Em. Max: 515-507/513-558 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

ZP-1 is supplied as a crystalline solid. A stock solution may be made by dissolving the ZP-1 in the solvent of choice. ZP-1 is soluble in the organic solvent dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of ZP-1 in DMF is approximately 0.5 mg/ml.

Description

Zinpyr-1 is a cell-permeable, fluorescein-based probe for zinc detection.¹ It undergoes a shift in excitation from 515 to 507 nm with increasing zinc concentrations using an emission filter over 513-558 nm.^{1,2}

References

1. Walkup, G.K., Burdette, S.C., Lippard, S.J., *et al.* A new cell-permeable fluorescent probe for Zn²⁺. *J. Am. Chem. Soc.* **122(23)**, 5644-5645 (2000).
2. Burdette, S.C., Walkup, G.K., Spingler, B., *et al.* Fluorescent sensors for Zn²⁺ based on a fluorescein platform: synthesis, properties and intracellular distribution. *J. Am. Chem. Soc.* **123(32)**, 7831-7841 (2001).

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, 10/25/2022

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