

PRODUCT INFORMATION



AlZin2

Item No. 40399

Formal Name: 5-((2-(2-((4-((bis(pyridin-2-ylmethyl)amino)methyl)-1H-imidazole-1-carbonyl)oxy)ethoxy)ethyl)carbamoyl)-3-oxo-3H-spiro[isobenzofuran-1,9'-xanthene]-3',6'-diyl diacetate

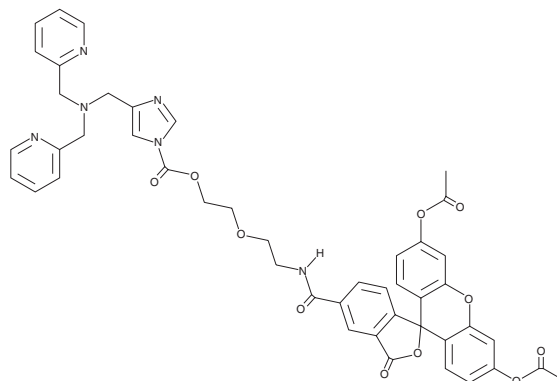
MF: C₄₆H₄₀N₆O₁₁

FW: 852.8

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

AlZin2 is supplied as a solid. A stock solution may be made by dissolving the AlZin2 in the solvent of choice, which should be purged with an inert gas. AlZin2 is soluble DMSO.

Description

AlZin2 is a conditional labeling compound for zinc-binding proteins that contains a zinc-coordinating tridentate dipicolylamine group, an electrophile-reactive acyl imidazole moiety, and fluorescein diacetate (Item No. 35701).¹ It is attached to proteins *via* an electrophilic attack on the acyl imidazole moiety when coordinated with protein-associated zinc. AlZin2 has been used to label zinc-binding proteins in HeLa cells for fluorescent detection and for LC-MS/MS proteomic experiments. It displays an emission at 559 nm upon excitation at 473 nm.

Reference

1. Miki, T., Awad, M., Nishikawa, Y., *et al.* A conditional proteomics approach to identify proteins involved in zinc homeostasis. *Nat. Methods* **13**(11), 931-937 (2016).

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, 07/01/2025

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