

ZnAF-2 . tetrahydrochloride

Fluorescent probe for Zn²⁺

Non cell permeable fluorescent reagent (Ex(max): 492nm; Em(max): 514nm) for the detection of low concentration of zinc ion due to its strong affinity to zinc ion (dissociation constant: 2.7nM). The sample zinc ion can be specifically detected. Low background fluorescence supersensitizes the visualization for *in vivo* sample zinc ion. Also available as cell permeable derivative (Prod. No. ALX-620-076). ZnAF-2 conjugates TPEN (Prod. No. ALX-450-011) affinities with fluorescein and enhances the specificity for Zn²⁺.

Zinc (Zn) is the second most abundant transition metal in the body and it is essential as catalytic, structural and regulatory ion. Zinc ions are involved in homeostasis, immune responses, oxidative stress, apoptosis and aging. Zinc has been proposed to function as a conventional neurotransmitter for the presynaptic neuron and as a transmembrane signal to traverse the postsynaptic neuron. Aberrant zinc metabolism is associated with many neurological diseases including Alzheimer's disease, Parkinson's disease and epilepsy. The most suitable technique for *in vivo* monitoring of zinc has been proven to be fluorescent imaging.

Citations: 3

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

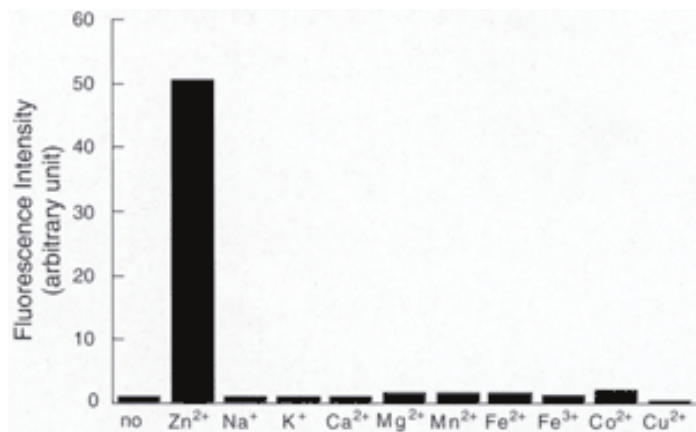


Figure 4: Changes in the fluorescence intensity of ZnAF-2 by adding each type of metal ion. 5 μ M of ZnAF-2 was dissolved in 100mM HEPES, pH 7.4. The fluorescence intensity when no metal ion was added was 1 (the concentration of the added metal ion was: 5 μ M Zn²⁺, Mn²⁺, Fe²⁺, Fe³⁺, Co²⁺ and Cu²⁺, 5mM of Na⁺, K⁺, Ca²⁺ and Mg²⁺)

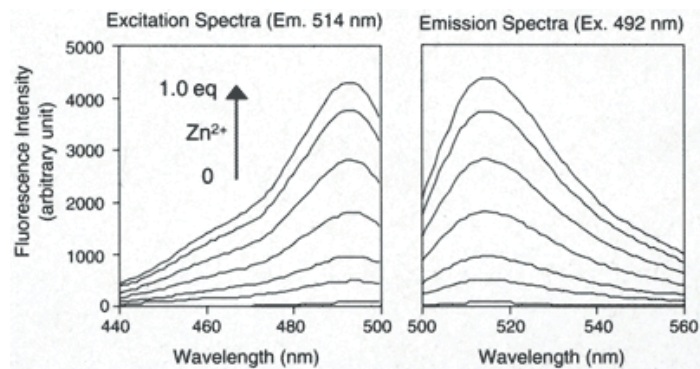


Figure 3: Excitation and fluorescence spectrum of ZnAF-2. 5 μ M of ZnAF-2 was dissolved in 100mM HEPES, pH 7.4. Each sample was 0, 0.5, 1, 2, 3, 4 and 5 μ M of zinc sulfate.

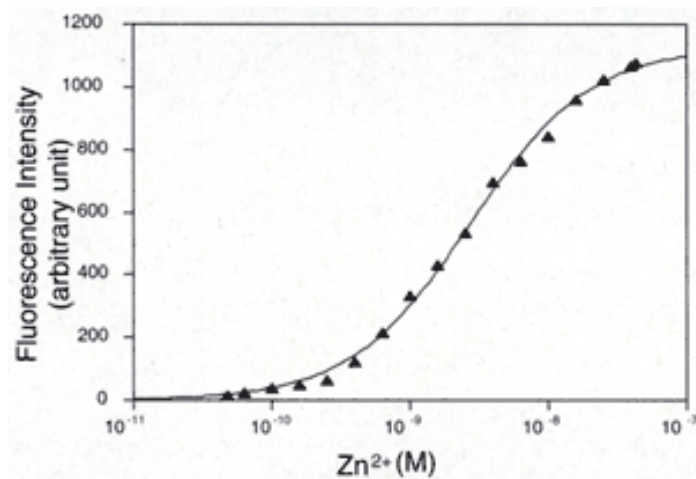


Figure 2: Changes in the fluorescence intensity of ZnAF-2 by adding Zn²⁺. 1 μ M of ZNAF-2 was dissolved in 100mM HEPES, pH 7.4.

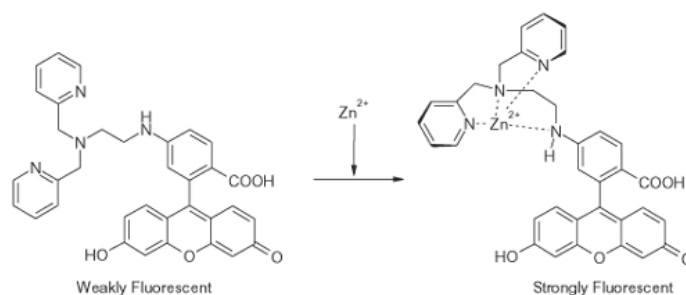
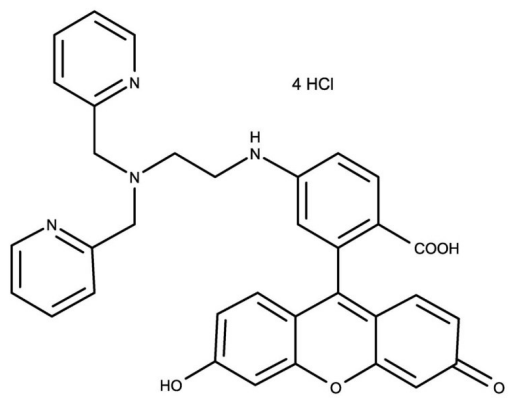


Figure 1: Reaction of ZnAF-2 with Zn²⁺.



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Prepare 500-5'000-fold dilution (~1-10µM) in 100mM HEPES, pH 7.4 immediately before use. Do not store diluted solutions. BSA and phenol red may affect the fluorescence and must be avoided.
Handling	After opening, prepare aliquots and store at +4°C. Protect from light. Keep under inert gas.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	6-[N-[N',N'-bis(2-Pyridinylmethyl)-2-aminoethyl]amino-3',6'-dihydroxy-spiro[isobenzofuran-1(3H),9'-[9H]xanthen]-3-one . 4 HCl
Appearance	Pale yellow liquid.
Concentration	~5mM
Formula	$C_{34}H_{28}N_4O_5 \cdot 4HCl$
Formulation	Dissolved in 0.28ml DMSO.
MW	572.6 . 145.8
Purity	≥98% (HPLC)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com