

## Fluorescent Indicators for Zinc

**Table 1.** Contents and storage information.

Material	Amount*	Storage	Stability
Water-soluble indicator salts	1 mg or 500 µg	<ul style="list-style-type: none"> <li>• ≤−20°C</li> <li>• Desiccate</li> <li>• Protect from light</li> </ul>	When stored as directed, this product is stable for at least 6 months.
Acetoxymethyl (AM) and acetate ester derivatives	1 mg, 100 µg, or 50 µg		
*See product label for details. <b>Note:</b> Avoid repeated freezing and thawing of DMSO stock solutions.			
<b>Approximate fluorescence excitation/emission maxima:</b> See Table 2.			

## Introduction

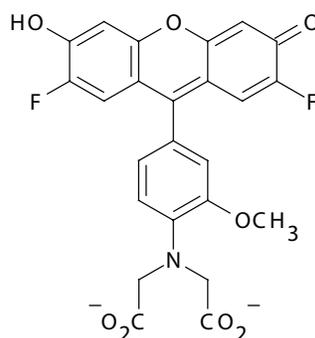
Zinc is an important divalent cation in biological systems, influencing DNA synthesis, microtubule polymerization, gene expression, apoptosis, immune system function, and the activity of enzymes such as carbonic anhydrase and matrix metalloproteinases (MMP). Zn<sup>2+</sup> is also functionally active in synaptic transmission, and is a contributory factor in neurological disorders including epilepsy and Alzheimer's disease.<sup>1,2</sup> The intracellular concentration of free Zn<sup>2+</sup> is extremely low in most cells (<1 nM), with the remainder being bound to proteins or nucleic acids.<sup>3</sup>

Zinc concentrations in the 1–100 nM range can be measured using fluorescent indicators nominally designed for Ca<sup>2+</sup> detection (*e.g.*, fura-2,  $K_d(\text{Zn}^{2+}) = 3 \text{ nM}$ ).<sup>4–6</sup> The FluoZin™-1 and FluoZin™-2 indicators are designed for detection of higher Zn<sup>2+</sup> concentrations that are present in synaptic vesicles and released in response to electrical stimulation or excitotoxic agonists.<sup>1,7</sup> These indicators based on the N-(2-methoxyphenyl)iminodiacetate chelator (Figure 1) are designed for detection of Zn<sup>2+</sup> in the 0.05–50 µM range with minimal interfering Ca<sup>2+</sup> sensitivity. FluoZin™-3 ( $K_d(\text{Zn}^{2+}) \sim 15 \text{ nM}$ ) is suitable for detection of Zn<sup>2+</sup> concentrations in the 1–100 nM range and it has shown to be the most Zn<sup>2+</sup>-sensitive and Zn<sup>2+</sup>-specific of the three FluoZin™ Zn<sup>2+</sup> indicators. The cell permeant AM-ester is useful for detecting low intracellular Zn<sup>2+</sup> levels and small concentration changes.<sup>19</sup>

RhodZin™-3 shows promise as an indicator for Zn<sup>2+</sup> in the mitochondria.<sup>20</sup> Exhibiting pH-insensitive fluorescence similar to tetramethylrhodamine, this indicator may prove useful for monitoring changes in Zn<sup>2+</sup> levels associated with pro-apoptosis.

The Newport Green™ DCF indicator has moderate zinc-binding affinity ( $K_d(\text{Zn}^{2+}) \sim 1 \text{ µM}$ ) but is essentially insensitive to Ca<sup>2+</sup> ( $K_d(\text{Ca}^{2+}) > 100 \text{ µM}$ ), making this a valuable probe for detecting Zn<sup>2+</sup> influx into neurons through voltage- or glutamate-gated channels.<sup>8–11</sup> When used alongside dyes with dual Ca<sup>2+</sup>/Zn<sup>2+</sup> sensitivity such as fura-2 and mag-fura-2, Newport Green™ DCF provides confirmation that changes in Zn<sup>2+</sup> levels, and not Ca<sup>2+</sup> or Mg<sup>2+</sup>, are

being detected.<sup>8,12</sup> Newport Green™ PDX incorporates the same di-(2-picolyl)amine chelator as Newport Green™ DCF but has a higher Zn<sup>2+</sup> dissociation constant (Table 1) and a larger Zn<sup>2+</sup>-free to Zn<sup>2+</sup>-saturated fluorescence intensity increase.



**Figure 1.** Structure of FluoZin™-1.

Table 2. Fluorescent indicators for zinc.

Indicator	Water-soluble form	Permeant ester form	Ex/Em*	K <sub>d</sub> (Zn <sup>2+</sup> )**
FluoZin™-1	F24180	F24181	495/515	8 μM
FluoZin™-2	—	F24189	495/525	2 μM
FluoZin™-3	F24194	F24195	494/516	15 nM <sup>§</sup>
RhodZin™-3	R36350	R36351	550/575	65 nM
Newport™ Green DCF	N7990	N7991	505/535	1 μM
Newport™ Green PDX	N24190	N24191	495/520	30 μM

\*Ex/Em: Fluorescence excitation and emission maxima in nm. \*\*Dissociation constant of indicator/Zn<sup>2+</sup> complex measured in 50 mM MOPS pH 7.0 at 22°C. <sup>§</sup>K<sub>d</sub> determined in 135 mM NaCl, 1.1 mM EGTA, 20 mM HEPES, pH 7.4, 0 to 10 μM free Zn<sup>2+</sup> at 22°C.

**Properties** Indicator spectroscopic properties and Zn<sup>2+</sup> binding affinities are summarized in Table 2. The other indicators listed in Table 2 all exhibit fluorescence intensity increases on binding Zn<sup>2+</sup> with no accompanying spectral shift. All of the indicators have negligible sensitivity to Ca<sup>2+</sup> concentrations up to at least 100 μM.

## Before Starting

### Preparing Stock Solutions and Storage

Prepare the stock solutions of water-soluble indicator salts in distilled water or aqueous buffers, and store frozen (≤−20°C) and **protected from light**. When stored properly, the solutions are stable for at least six months.

Acetoxymethyl (AM) and acetate ester derivatives are susceptible to hydrolysis and should be stored at  $\leq -20^{\circ}\text{C}$ , desiccated, and **protected from light**. When stored under these conditions, these compounds are stable for at least six months. Reconstitute ester derivatives in dimethylsulfoxide (DMSO). Concentrations of about 1–5 mM (molecular weights (MW) are printed on the product label) are generally suitable for these stock solutions. Once prepared, use DMSO stock solutions as soon as possible to avoid decomposition and a resulting loss of cell loading efficacy. Store stock solutions of AM esters frozen, desiccated, and **protected from light**. Because the integrity of AM esters is primarily dependent on minimizing their exposure to water, we recommend that you use high-quality anhydrous DMSO. **Avoid repeated freezing and thawing of DMSO stock solutions.**

## Experimental Protocols

---

### Cell Loading Guidelines

You may load the water-soluble salt forms of fluorescent indicators into cells by microinjection (see **Note** below) or by diffusion from a patch pipette.<sup>13</sup>

**Note:** Typical injection volume of 1–10 mM indicator solution is ~1% of the cell volume, giving a final intracellular indicator concentration of 10–100  $\mu\text{M}$ .

The following loading protocols using cell-permeant AM esters are provided as an introductory guide only; refer to published procedures for details.<sup>14–16</sup>

- 1.1 Dilute an aliquot of DMSO stock solution (1–5 mM) to a final concentration of 1–5  $\mu\text{M}$  in the buffered physiological medium of choice.

Addition of non-ionic detergent Pluronic® F-127 can assist in dispersing the non-polar AM ester in aqueous media. You can conveniently accomplish this by mixing the aliquot of AM ester stock solution in DMSO with an equal volume of 20% (w/v) Pluronic in DMSO (Cat. no. P3000MP) before diluting in loading medium, making the final Pluronic concentration about 0.02%. Invitrogen offers Pluronic® F-127 in 30 mL units of a sterile 10% (w/v) solution in water (Cat. no. P6866) and 2 g solid units (Cat. no. P6867).

- 1.2 Incubate the cells with the AM ester for 15 to 60 minutes at  $20^{\circ}\text{C}$  to  $37^{\circ}\text{C}$ .

You need to empirically determine the exact loading concentration, time, and temperature; in general it is desirable to use the minimum dye concentration required to yield fluorescence signals with adequate signal to noise levels. Subcellular compartmentalization, an inherent problem with the AM ester loading technique, is usually lessened by lowering the incubation temperature.<sup>14,15</sup>

- 1.3 Before you begin fluorescence measurements, wash cells in indicator-free medium to remove any dye that is nonspecifically associated with the cell surface, and then incubate for a further 30 minutes to allow complete de-esterification of intracellular AM esters.

### Response Calibration

You can carry out response calibration by measuring the fluorescence intensity of the carboxylate salt form of the indicator in solutions with precisely known free  $\text{Zn}^{2+}$  concentrations. If the  $\text{Zn}^{2+}$  concentrations are unbuffered, the approximation  $\text{Zn}^{2+}_{\text{total}} \sim \text{Zn}^{2+}_{\text{free}}$  is only valid if the indicator concentration is very low ( $< 0.1 \mu\text{M}$ ). You can prepare EGTA-buffered zinc calibration solutions using analogous calcium buffer methodology.<sup>17,18</sup>  $\text{Zn}^{2+}$  concentrations from 1 nM–10  $\mu\text{M}$  have been obtained in this way for calibration of the indicator mag-fura-5.<sup>5</sup> Chelators with lower  $\text{Zn}^{2+}$  affinity such as acetamidoiminodiacetic acid (ADA;  $K_d(\text{Zn}^{2+}) = 1.27 \times 10^{-7} \text{ M}$  at pH 7,  $I = 0.1$  and  $20^{\circ}\text{C}$ ) provide more optimal buffering in the micromolar range than EGTA ( $K_d(\text{Zn}^{2+}) = 6.8 \times 10^{-9} \text{ M}$  at pH 7,  $I = 0.1$  and  $20^{\circ}\text{C}$ ).

For indicators such as FluoZin™-1 and Newport Green™ DCF, you may determine the  $K_d$  using the following equation, in which F denotes fluorescence intensity measured at a single wavelength:

$$[Zn^{2+}] = K_d \frac{(F - F_{\min})}{(F_{\max} - F)}$$

In the above equation, the values of F are dependent on the concentration of indicator. It is important to recognize that the ion-binding and spectroscopic properties of fluorescent indicators can vary quite markedly in cellular environments. Consequently, *in situ* response calibrations of intracellular indicators often yield  $K_d$  values significantly different from *in vitro* determinations. Perform *in situ* calibrations of zinc indicators by exposing loaded cells to controlled  $Zn^{2+}$  concentrations in the presence of an ionophore. Pyrithione at a working concentration of 20  $\mu$ M is the most widely used  $Zn^{2+}$  ionophore.<sup>5,8,9</sup> Alternatively, 4-bromo A-23187 (Cat. no. B1494) has been used in some cases.<sup>6</sup> The zero reference level ( $F_{\min}$ ) for intracellular  $Zn^{2+}$  calibrations is usually set by adding 50–100  $\mu$ M TPEN (Cat. no. T-1210;  $K_d(Zn^{2+}) = 2.6 \times 10^{-16}$  M).

## References

1. Trends Pharmacol Sci 21, 395 (2000);
2. Science 265, 1464 (1994);
3. J Membrane Biol 123, 63 (1991);
4. J Biol Chem 270, 2473 (1995);
5. J Neurosci 17, 9554 (1997);
6. J Neurochem 71, 2401 (1998);
7. Nature 308, 734 (1984);
8. Proc Natl Acad Sci 96, 2414 (1999);
9. J Neurosci 19, RC31 (1999);
10. Eur J Neurosci 12, 3813 (2000);
11. J Physiol 528, 39 (2000);
12. J Neurochem 75, 1878 (2000);
13. *Imaging Neurons: A Laboratory Manual* R. Yuste, F. Lanni and A. Konnerth, Eds., pp 35.1–35.10, Cold Spring Harbor Laboratory Press (2000);
14. Methods Enzymol 302, 341 (1999);
15. Methods Enzymol 307, 441 (1999);
16. *Cell Biology: A Laboratory Handbook, 2<sup>nd</sup> Edition*, J.E. Celis, Ed., Volume 3, pp 363–374, Academic Press (1998);
17. Methods Cell Biol 40, 3 (1994);
18. Methods Enzymol 172, 230 (1989);
19. Cell Calcium 31, 245 (2002);
20. Cell Calcium 34, 281 (2003).

## Product List Current prices may be obtained from our website or from our Customer Service Department.

Cat. no.	Product Name	Unit Size
B1494	4-bromo A-23187, free acid .....	1 mg
F24180	FluoZin™-1, tripotassium salt *cell impermeant* .....	500 $\mu$ g
F24181	FluoZin™-1, AM *cell permeant* .....	50 $\mu$ g
F24189	FluoZin™-2, AM *cell permeant* .....	50 $\mu$ g
F24194	FluoZin™-3, tetrapotassium salt *cell impermeant* .....	500 $\mu$ g
F24195	FluoZin™-3, AM *cell permeant* .....	100 $\mu$ g
N24190	Newport Green™ PDX .....	1 mg
N24191	Newport Green™ PDX acetoxymethyl ether.....	1 mg
N7990	Newport Green™ DCF, dipotassium salt *cell impermeant* .....	1 mg
N7991	Newport Green™ DCF diacetate *cell permeant* .....	1 mg
P10020	PowerLoad™ concentrate, 100X .....	5 mL
P3000MP	Pluronic® F-127 *20% solution in DMSO* .....	1 mL
P6866	Pluronic® F-127 *10% solution in water* *0.2 $\mu$ m filtered* .....	30 mL
P6867	Pluronic® F-127 *low UV absorbance* .....	2 g
R36350	RhodZin™-3, dipotassium salt *cell impermeant* .....	500 $\mu$ g
R36351	RhodZin™-3, AM *cell permeant* .....	50 $\mu$ g
T1210	tetrakis-(2-pyridylmethyl)ethylenediamine (TPEN) .....	100 mg

## Contact Information

---

### **Molecular Probes, Inc.**

29851 Willow Creek Road  
Eugene, OR 97402  
Phone: (541) 465-8300  
Fax: (541) 335-0504

### **Customer Service:**

6:00 am to 4:30 pm (Pacific Time)  
Phone: (541) 335-0338  
Fax: (541) 335-0305  
probesorder@invitrogen.com

### **Toll-Free Ordering for USA:**

Order Phone: (800) 438-2209  
Order Fax: (800) 438-0228

### **Technical Service:**

8:00 am to 4:00 pm (Pacific Time)  
Phone: (541) 335-0353  
Toll-Free (800) 438-2209  
Fax: (541) 335-0238  
probestech@invitrogen.com

### **Invitrogen European Headquarters**

Invitrogen, Ltd.  
3 Fountain Drive  
Inchinnan Business Park  
Paisley PA4 9RF, UK  
Phone: +44 (0) 141 814 6100  
Fax: +44 (0) 141 814 6260  
Email: euroinfo@invitrogen.com  
Technical Services: eurotech@invitrogen.com

**For country-specific contact information,  
visit [www.invitrogen.com](http://www.invitrogen.com).**

Further information on Molecular Probes products, including product bibliographies, is available from your local distributor or directly from Molecular Probes. Customers in Europe, Africa and the Middle East should contact our office in Paisley, United Kingdom. All others should contact our Technical Service Department in Eugene, Oregon.

Molecular Probes products are high-quality reagents and materials intended for research purposes only. These products must be used by, or directly under the supervision of, a technically qualified individual experienced in handling potentially hazardous chemicals. Please read the Material Safety Data Sheet provided for each product; other regulatory considerations may apply.

### **Limited Use Label License No. 223: Labeling and Detection Technology**

The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) to not transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. Invitrogen Corporation will not assert a claim against the buyer of infringement of the above patents based upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Invitrogen is willing to accept return of the product with a full refund. For information on purchasing a license to this product for purposes other than research, contact Molecular Probes, Inc., Business Development, 29851 Willow Creek Road, Eugene, OR 97402, Tel: (541) 465-8300. Fax: (541) 335-0354.

Several Molecular Probes products and product applications are covered by U.S. and foreign patents and patents pending. All names containing the designation ® are registered with the U.S. Patent and Trademark Office.

Copyright 2009, Molecular Probes, Inc. All rights reserved. This information is subject to change without notice.