

Metal Fluor™ Zn-520, Potassium Salt

 Catalog number: 21262
 Unit size: 1 mg

Component	Storage	Amount (Cat No. 21262)
Metal Fluor™ Zn-520, Potassium Salt	Freeze (< -15 °C), Minimize light exposure	1 vial (1 mg)

OVERVIEW

Metal Fluor™ Zn-520 is designed for detection of higher zinc ion concentrations that are present in synaptic vesicles and released in response to electrical stimulation or excitotoxic agonists (0.05-50 μM) with minimal interfering calcium sensitivity. Metal Fluor™ Zn-520 has shown great fluorescence enhancement upon binding zinc ion (>250 folds). This cell-impermeant zinc ion indicator is useful for detecting extracellular zinc ion.

AT A GLANCE
Protocol Summary

1. Test samples (50 μL) or Zn²⁺ Standard
2. Add Zinc Probe Reagent (50 μL)
3. Incubate at room temperature for 5 - 10 minutes
4. Read fluorescence at Ex/Em= 485/525 nm

Important Note

Thaw the dye at room temperature before starting the experiment.

KEY PARAMETERS
Fluorescence microplate reader

Cutoff	515 nm
Emission	525 nm
Excitation	485 nm
Recommended plate	Solid black

PREPARATION OF STOCK SOLUTIONS

Unless otherwise noted, all unused stock solutions should be divided into single-use aliquots and stored at -20 °C after preparation. Avoid repeated freeze-thaw cycles

Zinc standard stock solution (1 mM)

Add 10 μL of 100 mM ZnCl₂ Standard into 990 μL HEPES buffer to get 1 mM ZnCl₂ standard solution.

Metal Fluor™ Zn-520

Make 1-4 mM stock solution in water.

Note: Store at -20°C in single-use aliquots.

PREPARATION OF STANDARD SOLUTIONS

For convenience, use the Serial Dilution Planner:
<https://www.aatbio.com/tools/serial-dilution/21262>

ZnCl₂ standard

Add 100 μL of 1 mM Zinc standard solution to 900 μL HEPES buffer to get 100 μM ZnCl₂ standard solution (Zn7). Then take the 100 μM ZnCl₂ standard solution to perform 1:3 serial dilutions to get serially diluted ZnCl₂ standards (Zn6 - Zn1).

PREPARATION OF WORKING SOLUTION
Metal Fluor™ Zn-520 working solution

Make a 10-50 μM working solution from the Metal Fluor™ Zn-520 stock solution.

Note: Working concentration may be optimized empirically.

SAMPLE EXPERIMENTAL PROTOCOL

Table 1. Layout of ZnCl₂ standards and test samples in a solid black 96-well microplate. Zn = Zinc standard (Zn1 - Zn7, 0.1 to 100 μM); BL = blank control; TS = test sample.

BL	BL	TS	TS
Zn1	Zn1
Zn2	Zn2
Zn3	Zn3		
Zn4	Zn4		
Zn5	Zn5		
Zn6	Zn6		
Zn7	Zn7		

Table 2. Reagent composition for each well.

Well	Volume	Reagent
Zn1 - Zn7	50 μL	Serial dilution (0.1 to 100 μM)
BL	50 μL	HEPES buffer
TS	50 μL	Sample

1. Dilute the test sample to a 5 - 100 μM range with HEPES buffer.
2. Prepare ZnCl₂ standards (Zn), blank controls (BL), and test samples (TS) according to the layout provided in Table 1 and Table 2. For a 384-well plate, use 25 μL of reagent per well instead of 50 μL.
3. Add 50 μL of Zinc working solution into each well of ZnCl₂ standard, blank control, and test samples to make the total ZnCl₂ assay volume of 100 μL/well. For a 384-well plate, add 25 μL of HEPES buffer into each well instead, for a total volume of 50 μL/well.
4. Incubate the reaction for 5 - 10 minutes at room temperature, protected from light.
5. Monitor the fluorescence increase with a fluorescence plate reader at Ex/Em = 485/525 nm.

EXAMPLE DATA ANALYSIS AND FIGURES

The reading (RFU) obtained from the blank standard well is used as a negative control. Subtract this value from the other standards' readings to obtain the baseline corrected values. Then, plot the standards' readings to obtain a standard curve and equation. This equation can be used to calculate Zinc Chloride samples. We recommend using the Online Linear Regression Calculator which can

be found at:

<https://www.aatbio.com/tools/linear-logarithmic-semi-log-regression-online-calculator>

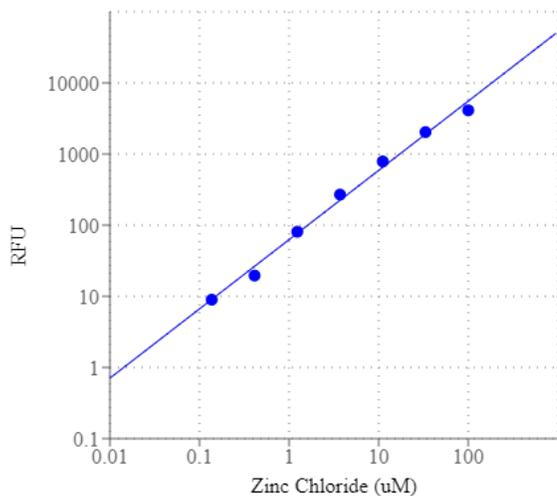


Figure 1. Zinc Chloride dose response was measured on a 96-well black plate with the Metal Fluor™ Zn-520.

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.