

FDP [Fluorescein diphosphate, tetraammonium salt] *CAS 217305-49-2*

 Catalog number: 11600, 11601
 Unit size: 5 mg, 10 mg

Component	Storage	Amount (Cat No. 11600)	Amount (Cat No. 11601)
FDP [Fluorescein diphosphate, tetraammonium salt] *CAS 217305-49-2*	Freeze (< -15 °C), Minimize light exposure	1 vial (5 mg)	1 vial (10 mg)

OVERVIEW

Upon interaction with phosphatases the colorless and non-fluorescent FDP is hydrolyzed to highly fluorescent fluorescein, which exhibits excellent spectral properties that match the optimal detection window of most fluorescence instruments that are equipped with the Argon laser excitation. Alternatively, FDP can also be used to detect phosphatases in a chromogenic mode since the enzymatic product (fluorescein) exhibits a large extinction coefficient (close to 100,000 cm⁻¹mol⁻¹). In some literature, FDP was considered to be one of the most sensitive fluorogenic phosphatase substrates. FDP has been widely used in various ELISA assays. Additionally it is also used to detect tyrosine phosphatases. FDP is thermally unstable, and special cautions need be excised for storing the solid sample and stock solutions.

AT A GLANCE
Protocol Summary

1. Prepare 10 - 50 μM Phosphate in Tris buffer (50 μL)
2. Add phosphatase standards and/or test samples
3. Incubate at room temperature or 37°C for 30 to 120 minutes
4. Monitor fluorescence intensity at Ex/Em= 490/514 nm

Important Note

Phosphatase substrates can be detected by phosphatases that may not be specifically listed. The following is the recommended protocol for phosphatase assay in solution. The protocol only provides a guideline, should be modified according to the specific needs.

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

FDP stock solution

Prepare a 2 to 10 mM stock solution in ddH₂O.

Note: The stock solution should be used promptly.

PREPARATION OF WORKING SOLUTION
FDP working solution (2X)

On the day of the experiment, either dissolve FDP in ddH₂O or thaw an aliquot of the stock solution at room temperature. Prepare a 2X working solution of 10 to 50 μM in 100 mM Tris buffer or buffer of your choice, pH 8 to 9 (not phosphate buffer).

SAMPLE EXPERIMENTAL PROTOCOL

1. Add 50 μL of 2X FDP working solution into each well of the phosphatase standard, blank control, and test samples to make the total phosphatase assay volume of 100 μL/well. For a 384-well plate, add 25 μL of sample and 25 μL of 2X Phosphate working solution into each well.
2. Incubate the reaction for 30 to 120 minutes at the desired

temperature, protected from light.

3. Monitor the fluorescence increase at an appropriate filter set with a fluorescence plate reader.
4. The fluorescence in blank wells (with the assay buffer only) is used as a control, and is subtracted from the values for those wells with the phosphatase reactions.

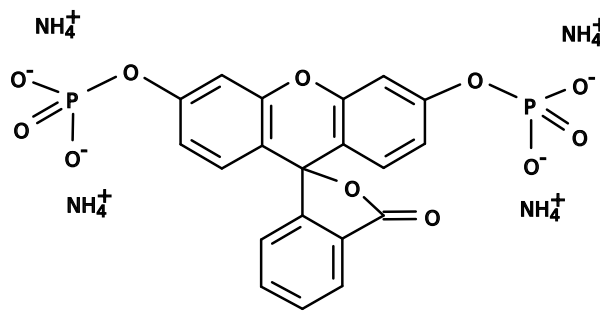
EXAMPLE DATA ANALYSIS AND FIGURES


Figure 1. Chemical structure for FDP [Fluorescein diphosphate, tetraammonium salt] *CAS 217305-49-2*

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.