

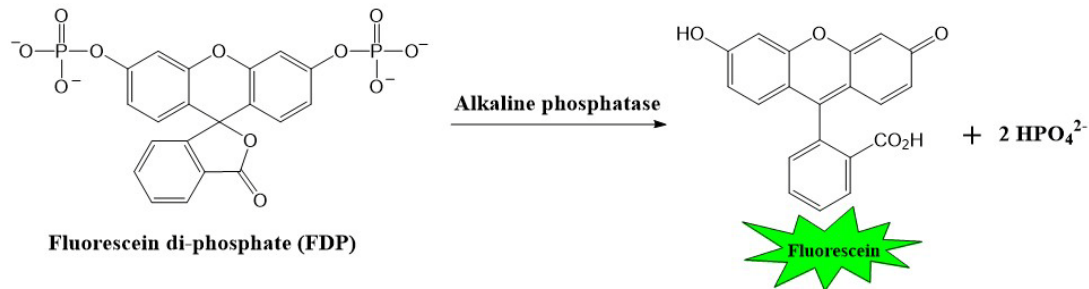
FDP (Fluorescein diphosphate, tetraammonium salt)

Catalog Number	Packaging Size
C275	2 mg

Storage upon receipt: -20°C, protected from light

Introduction

FDP is a colorless and nonfluorescent substrate for alkaline phosphatases. Sequential alkaline phosphatase mediated hydrolysis of its two phosphate substituents yields weakly fluorescent fluorescein monophosphate followed by strongly fluorescent fluorescein (excitation/emission ~488/515 nm).



Specifications

Label:	Fluorescein	
Ex/Em:	488/515 nm	
Detection Method:	Fluorescent	
Molecular Formula:	$\text{C}_{20}\text{H}_{26}\text{N}_4\text{O}_{11}\text{P}_2$	
Molecular Weight:	560.39	
CAS Number:	217305-49-2	
Storage Conditions:	-20°C, protect from light	
Shipping Condition:	Room Temperature	

Applications

Phosphatase Substrate

References:

1. Evaluation of fluorescent compound interference in 4 fluorescence polarization assays: 2 kinases, 1 protease, and 1 phosphatase.

- Turek-Etienne TC, Small EC, Soh SC, Xin TA, Gaitonde PV, Barrabee EB, Hart RF, Bryant RW
J Biomol Screen (2003) 8:176-176
2. Fluorogenic substrates for beta-D-galactosidases and phosphatases derived from fluorescein (3,6-dihydroxyfluoran) and its monomethylether. Fluorogenic substrates for beta-D-galactosidases and phosphatases derived from fluorescein (3,6-dihydroxyfluoran) and its monomethylether.
ROTMAN B, ZDERIC JA, EDELSTEIN M,
Proc Natl Acad Sci U S A (1963) 50:1-6
 3. Ionization and tautomerism of oxyxanthene dyes in aqueous butanol ionization and tautomerism of oxyxanthene dyes in aqueous butanol
Mchedlov-Petrosyan NO, Tychina ON, Berezhnaya TA, Alekseeva VI, Savvina LP
Dyes Pigment (1999) 43:33-46