

FDP (high purity)

FDP (fluorescein diphosphate, tetraammonium salt, high purity) is the most sensitive fluorogenic alkaline phosphatase substrate. It is also an excellent substrate for protein tyrosine phosphatase (PTPase) and thus the substrate of choice in the high throughput screening (HTS) for PTPase inhibitor.



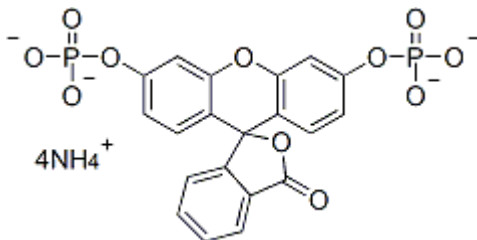
Product attributes

CAS number	217305-49-2
Excitation/Emission	490/514 nm (end product)

Product Description

FDP (fluorescein diphosphate, tetraammonium salt, high purity) is the most sensitive fluorogenic alkaline phosphatase substrate. It is also an excellent substrate for protein tyrosine phosphatase (PTPase) and thus the substrate of choice in the high throughput screening (HTS) for PTPase inhibitor. FDP itself is colorless, but the end product of the enzymatic reaction is the green fluorescent dye fluorescein.

- $\lambda_{Ex}/\lambda_{Em} = 490/514$ nm (end product)
- $\epsilon = 90,000$ (end product)
- Off-white solid soluble in water
- Store at -20°C and protect from light, especially in solution
- $\text{C}_{20}\text{H}_{26}\text{N}_4\text{O}_{11}\text{P}_2$
- MW: 560.39
- [217305-49-2]



References

1. Appl Environ Microbiol 64, 1526 (1998).
2. Biol Res 27, 241 (1994).
3. Anal Biochem 285, 58 (2000).
4. J Biomol Screen 4, 327 (1999).
5. Proc Natl Acad Sci USA 93, 3068 (1996).

This datasheet was generated on June 5, 2026 at 02:37:36 AM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/fdp-fluorescein-diphosphate-tetraammonium-salt-high-purity/>