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.

alictum &

Product attributes

CAS number 217305-49-2

Excitation/Emission 490/514 nm (end product)

Call us: 800-304-5357 Email: btinfo@biotium.com

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)
- $\varepsilon = 90,000$ (end product)
- Off-white solid soluble in water
- Store at -20°C and protect from light, especially in solution
- C20H26N4O11P2
- 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 November 4, 2025 at 06:47:36 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/fdp-fluorescein-diphosphate-tetraammonium-salt-high-purity/