

## MUP, Na

MUP, Na (4-Methylumbelliferyl phosphate, disodium salt) is the sodium salt of MUP. MUP sodium salt and MUP free acid (#10009) are the most widely used fluorogenic substrates for alkaline phosphatase detection.



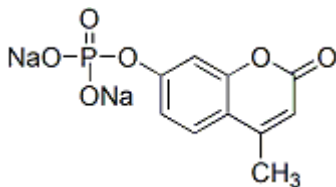
### Product attributes

CAS number	22919-26-2
Excitation/Emission	360/450 nm (end product)

## Product Description

MUP, Na (4-Methylumbelliferyl phosphate, disodium salt) is the sodium salt of MUP. MUP sodium salt and MUP free acid (#[10009](#)) are the most widely used fluorogenic substrates for alkaline phosphatase detection.

- $C_{10}H_7Na_2O_6P \cdot 3H_2O$
- MW: 354.2
- [22919-26-2]



This datasheet was generated on June 30, 2026 at 02:55:31 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/mupna-4-methylumbelliferyl-phosphate-disodium-salt-2/>