



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

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.



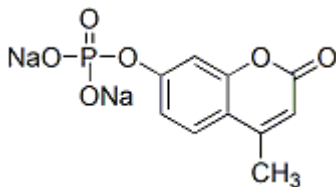
Product attributes

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 January 2, 2026 at 04:54:38 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/mupna-4-methylumbelliferyl-phosphate-disodium-salt-2/>