

Aminoxy-5(6)-FAM

Our aminoxy reagents such as aminoxy-5(6)-FAM react with aldehydes or ketones to form a stable oxime linkage under mild conditions, and can be used to label abasic sites in damaged DNA.



Product attributes

Excitation/Emission	494/521 nm (pH 9)
---------------------	-------------------

Product Description

Our aminoxy reagents such as aminoxy-5(6)-FAM react with aldehydes or ketones to form a stable oxime linkage under mild conditions. Similar to ARP (#90073), aminoxy-5(6)-FAM can be used to label abasic sites in damaged DNA. It can also be used to label polysaccharides and glycoproteins.

- $\lambda_{Ex}/\lambda_{Em}$ (pH 9) = 494/521 nm
- Orange solid soluble in DMF, DMSO, or pH >6.5 buffers
- Store at -20 °C and protect from moisture and light
- $C_{26}H_{21}F_3N_2O_9$
- MW: 562.5

