

## Aminoxy-5(6)-TAMRA

Aminoxy-5(6)-TAMRA reacts with aldehydes or ketones to form a stable oxime linkage under mild conditions. It can also be used to label abasic sites in damaged DNA.



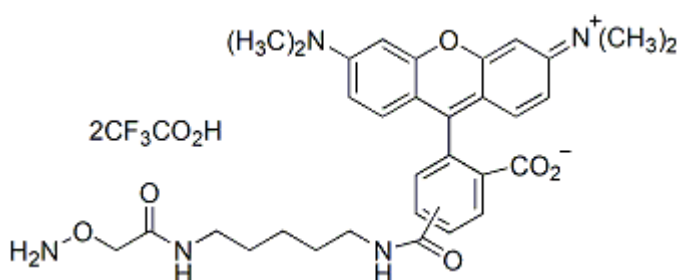
### Product attributes

Excitation/Emission	540/565 nm
Chemical reactivity (reacts with)	Aldehydes/ketones
Functional group	Aminoxy (hydroxylamine)
Storage Conditions	Store at -10 to -35 °C, Protect from light

## Product Description

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

- $\lambda_{Ex}/\lambda_{Em}(\text{MeOH}) = 540/565 \text{ nm}$
- Dark red solid soluble in DMF or DMSO
- Store at -20 °C and protect from moisture and light
- $\text{C}_{36}\text{H}_{39}\text{F}_6\text{N}_5\text{O}_{10}$
- MW: 815.71



## References

1. Polym Chem. (2015) 6(31): 5683–5692. [DOI: 10.1039/C5PY00282F](https://doi.org/10.1039/C5PY00282F)

This datasheet was generated on December 5, 2025 at 02:12:02 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/aminoxy-56-tamra/>