

Aminoxy-5(6)-ROX

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



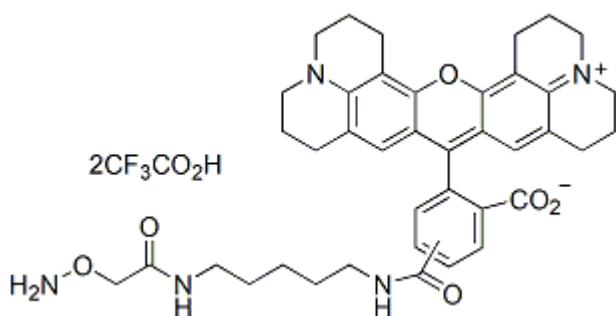
Product attributes

Excitation/Emission	568/595 nm
---------------------	------------

Product Description

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

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH): 568/595 nm
- Dark purple solid soluble in DMF or DMSO
- Store at -20 °C and protect from moisture and light
- C₄₅H₅₁F₆N₅O₁₀
- MW: 919.86



This datasheet was generated on June 3, 2026 at 09:30:58 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/aminoxy-6-rox/>