Cyanine Dye Aminooxy

Cyanine dyes with an aminooxy reactive group are useful for fluorescently labeling aldehyde or ketone groups on target molecules such as polysaccharids, glycoproteins or antibodies.



Product attributes

Chemical reactivity (reacts with)	Aldehydes/ketones
Functional group	Aminooxy (hydroxylamine)
Storage Conditions	Store at -10 to -35 °C. Protect from light

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

Cyanine dye with an aminooxy reactive group can be used to fluorescently label target molecules with an aldehyde or ketone group. Aminooxy group readily reacts with an aldehyde or ketone group to form a stable oxime linkage without the use of reducing agents. We offer Cyanine 647 Aminooxy and Cyanine 555 Aminooxy.

• React with aldehyde/ketone groups on target molecules.

Cyanine 555:

- λ_{Ex}/λ_{Em} (water) = 555/565 nm
- Dark blue solid soluble in water, DMF and DMSO
- C₃₈H₄₈F₆N₄O₁₂S₂
- MW: 931

$$HO_3S$$
 H
 $C=C-C$
 H
 N^+
 H
 $N^ O-NH_3^+$
 $2CF_3CO_2^-$

Cyanine 647:

- λ_{Ex}/λ_{Em} (water) = 650/665 nm
- Dark blue solid soluble in water, DMF and DMSO
- \bullet C₄₀H₅₀F₆N₄O₁₂S₂
- MW: 957

Also see our CF® Dye Aminooxy, available in a wide selection of our bright and photostable CF® Dyes.

This datasheet was generated on November 5, 2025 at 02:35:07 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/cyanine-dye-aminooxy/