

## Cyanine Dye Aminoxy

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



### Product attributes

<b>Chemical reactivity (reacts with)</b>	Aldehydes/ketones
<b>Functional group</b>	Aminoxy (hydroxylamine)
<b>Storage Conditions</b>	Store at -10 to -35 °C, Protect from light

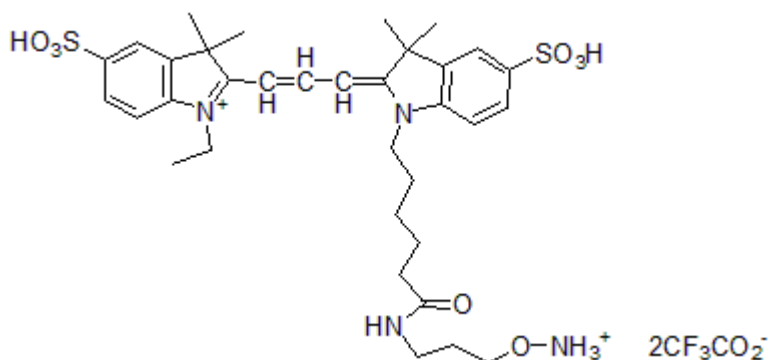
## Product Description

Cyanine dye with an aminoxy reactive group can be used to fluorescently label target molecules with an aldehyde or ketone group. Aminoxy 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 Aminoxy and Cyanine 555 Aminoxy.

- React with aldehyde/ketone groups on target molecules.

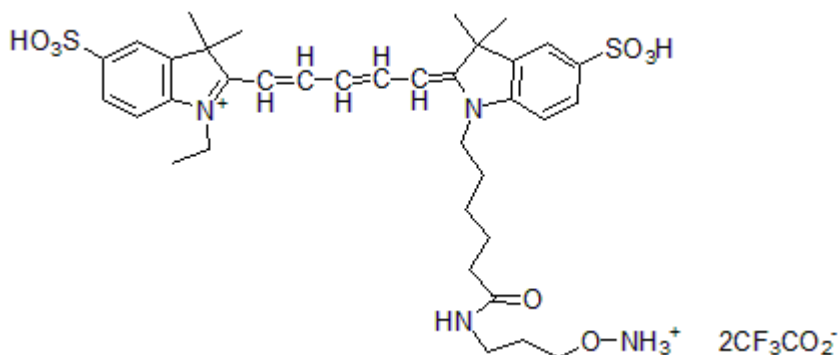
Cyanine 555:

- $\lambda_{Ex}/\lambda_{Em}$  (water) = 555/565 nm
- Dark blue solid soluble in water, DMF and DMSO
- $C_{38}H_{48}F_6N_4O_{12}S_2$
- MW: 931



Cyanine 647:

- $\lambda_{Ex}/\lambda_{Em}$  (water) = 650/665 nm
- Dark blue solid soluble in water, DMF and DMSO
- $C_{40}H_{50}F_6N_4O_{12}S_2$
- MW: 957



Also see our [CF® Dye Aminoxy](#), available in a wide selection of our bright and photostable CF® Dyes.