

Flubida-2

Flubida-2 has been used to detect pH at a specific site in a cell such as cell organelles by directing the probe to where avidin fusion proteins are located. The probe is a conjugate of biotin and fluorescein diacetate, which is nonfluorescent until the probe has entered the cells and hydrolyzed by endoesterases.



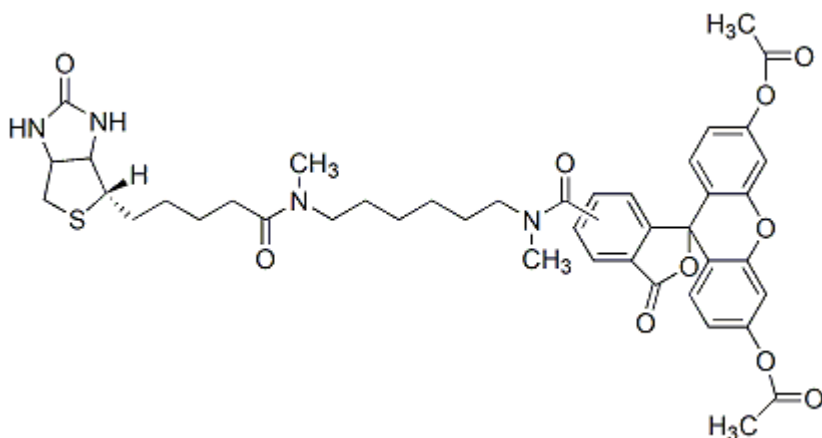
Product attributes

Cell permeability	Membrane permeant
pKa	pKa 6.5
Indicator type	Ratiometric
Colors	Green
Excitation/Emission	492/517 nm (after hydrolysis)

Product Description

Flubida-2 has been used to detect pH at a specific site in a cell such as cell organelles by directing the probe to where avidin fusion proteins are located. The probe is a conjugate of biotin and fluorescein diacetate, which is nonfluorescent until the probe has entered the cells and hydrolyzed by endoesterases. Flubida-2 is membrane-permeable and thus can be delivered into cells via simple incubation with the probe in a buffer.

- $\lambda_{Ex}/\lambda_{Em}$: 492/517 nm (pH 9) (after hydrolysis)
- Off-white solid soluble in DMSO
- Store at 4 °C
- $C_{43}H_{48}N_4O_9S$
- MW: 812.93



References

1. Chemistry and Biology 7, 197 (2000).

This datasheet was generated on January 19, 2026 at 11:14:07 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/flubida-2/>