

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

Call us: 800-304-5357

Cell permeability	Membrane permeant
рКа	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
- C₄₃H₄₈N₄O₉S
- MW: 812.93

References

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

This datasheet was generated on November 21, 2025 at 12:07:25 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/flubida-2/