Biotin-AM3-25

Biotin-AM3-25 is a conjugate of AM3-25 and biotin. Membranes labeled with the probe can subsequently couple to avidin or streptavidin.



Product attributes

Call us: 800-304-5357

	For live or fixed cells	For live/intact cells
	Fixation options	Fix after staining (formaldehyde), Permeab
	Conjugation	Biotin

Product Description

Biotin-AM3-25 is a conjugate of AM3-25 and biotin. Membranes labeled with the probe can subsequently couple to avidin or streptavidin.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 510/625 nm; $\lambda_{Ex}\lambda_{Em}$ (liposomes) = 480/600 nm
- Dark red soluble in MeOH or DMSO
- Store at -20°C and protect from light
- C₆₇H₁₁₈Cl₂N₆O₂S
- MW: 1142.66

$$[H_3C(H_2C)_{17}]_2N - CH = CH - CH_2)_3 - N_+(CH_2)_3NH - C - (CH_2)_4 - N_+(CH_2)_4 - N_+(CH_2)_4$$

This datasheet was generated on December 2, 2025 at 05:52:25 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/biotin-am3-25/