

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.



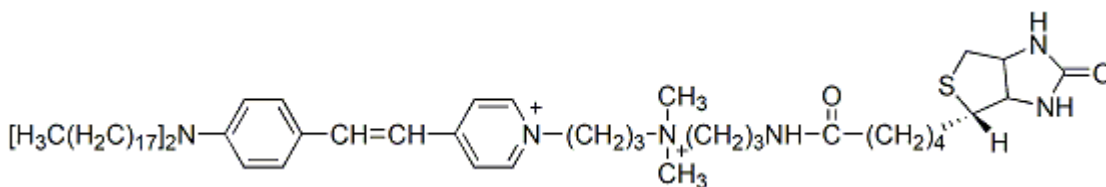
Product attributes

For live or fixed cells	For live/intact cells
Fixation options	Fix after staining (formaldehyde), Permeabilization
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
- $C_{67}H_{118}Cl_2N_6O_2S$
- MW: 1142.66



This datasheet was generated on January 24, 2026 at 01:10:09 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/biotin-am3-25/>