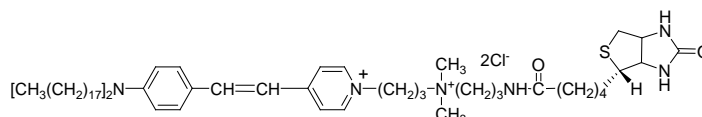


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Biotin AM3-25**CATALOG #:** 60032**MOLECULAR INFORMATION:** C₆₇H₁₁₈Cl₂N₆O₂S
Mwt: 1142.66**PROPERTIES:**

Color & Form	Dark red solid
Purity	≥ 95% by TLC
Solubility	Soluble in MeOH or DMSO
$\lambda_{ex} \backslash \lambda_{em}$	510 nm/ 625 nm (MeOH), 480 nm/ 600 nm (liposomes)
Extinction Coefficient	45,000 ± 3,000

STORAGE AND HANDLING: Store at -20 °C and protect from light.**APPLICATION:** Biotin-AM3-25 is a conjugate of AM3-25 and biotin. Membranes labeled with the probe can subsequently couple to avidin or streptavidin.**TOXICITY:** Unknown.

FIRST AID:	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.
-------------------	---

Disclaimer: <i>Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.</i>
--