Biotin-X-C5-maleimide

Biotin-X- C_5 -maleimide (N-(biotinoyl)-N'-(iodoacetyl)ethylenediamine) is readily and selectively coupled to thiols.



Product attributes

Call us: 800-304-5357

Product Description

 $Biotin-X-C_5-male imide \ (N-(biotinoyl)-N'-(iodoacetyl) ethylene diamine) \ is \ readily \ and \ selectively \ coupled \ to \ thiols.$

- White solid soluble in DMSO, or DMF
- Store at -20°C
- C₂₇H₃₉N₅O₇S
- MW: 521.67

This datasheet was generated on November 28, 2025 at 05:58:58 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/biotin-x-c5-maleimide-n-5-6-biotinoylaminohexanoylaminopentylmaleimide/