

## Biotin-X-C5-maleimide

Biotin-X-C5-maleimide (N-(biotinoyl)-N'-(iodoacetyl)ethylenediamine) is readily and selectively coupled to thiols.

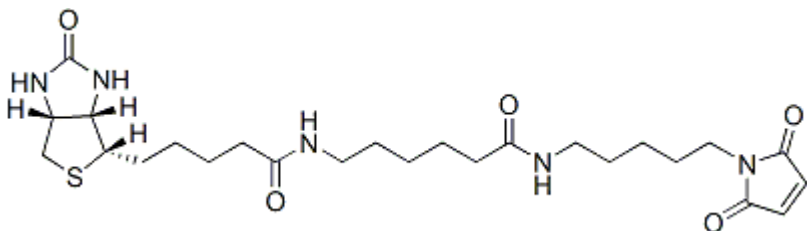


### Product attributes

## Product Description

Biotin-X-C5-maleimide (N-(biotinoyl)-N'-(iodoacetyl)ethylenediamine) is readily and selectively coupled to thiols.

- White solid soluble in DMSO, or DMF
- Store at -20 °C
- C<sub>27</sub>H<sub>39</sub>N<sub>5</sub>O<sub>7</sub>S
- MW: 521.67



This datasheet was generated on June 22, 2026 at 06:25:02 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/biotin-x-c5-maleimide-n-5-6-biotinoylamino-hexanoylamino-pentyl-maleimide/>