## **Biotin SE**

Biotin SE (Biotin, succinimidyl ester) is readily coupled to a primary or secondary amine under mild conditions.



#### **Product attributes**

Call us: 800-304-5357 Email: btinfo@biotium.com

# **Product Description**

Biotin SE (Biotin, succinimidyl ester) is readily coupled to a primary or secondary amine under mild conditions.

- White solid soluble in DMSO and moderately soluble in DMF
- Store at -20°C
- C<sub>14</sub>H<sub>19</sub>N<sub>3</sub>O<sub>5</sub>S
- MW: 341.38

### References

- 1. Bioconjugate Chem 8, 552 (1997).
- 2. Bioconjugate Chem 8, 545 (1997).
- 3. J Immunol 159, 5871 (1997).

This datasheet was generated on November 3, 2025 at 11:20:04 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/biotinsuccinimidyl-ester-biotin-se/">https://biotium.com/product/biotinsuccinimidyl-ester-biotin-se/</a>