## N-(1,10-Phenanthrolin-5-yl)iodoacetamide

N-(1,10-Phenanthrolin-5-yl)iodoacetamide has been used to label thiols to confer the metal-binding properties of the phenanthroline chelator. For example, completing of the chelator with copper can be used to selectively cleave RNA or DNA in the presence of hydrogen peroxide.



## **Product attributes**

Call us: 800-304-5357

## **Product Description**

N-(1,10-Phenanthrolin-5-yl)iodoacetamide has been used to label thiols to confer the metal-binding properties of the phenanthroline chelator. For example, completing of the chelator with copper can be used to selectively cleave RNA or DNA in the presence of hydrogen peroxide.

- Off-white solid soluble in DMF
- Store at -20°C
- C<sub>14</sub>H<sub>10</sub>IN<sub>3</sub>O
- MW: 363.15

## References

- 1. Nucleic Acids Res 22, 4789 (1994).
- 2. Bioconjugate Chem 4, 69 (1993).
- 3. Meth Enzymol 208, 414 (1991).
- 4. Anal Rev Biochem 59, 207 (1990).
- 5. J Am Chem Soc 111, 4941 (1989).
- 6. PNAS 86, 9702 (1989).

This datasheet was generated on November 30, 2025 at 12:06:12 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/n-110-phenanthrolin-5-yliodoacetamide/">https://biotium.com/product/n-110-phenanthrolin-5-yliodoacetamide/</a>