

## 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.

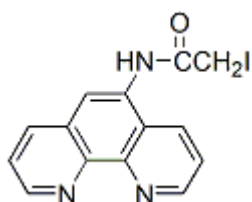


### Product attributes

### 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 January 20, 2026 at 03:06:55 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/n-110-phenanthrolin-5-yl-iodoacetamide/>