

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.

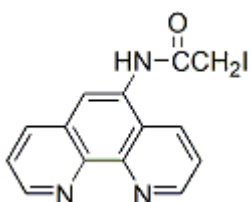


Product attributes

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
- C₁₄H₁₀IN₃O
- 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).