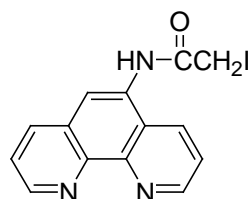


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: *N*-(1,10-Phenanthrolin-5-yl)iodoacetamide**CATALOG #** 92015**MOLECULAR INFORMATION:** C₁₄H₁₀IN₃O
MWt: 363.15**PROPERTIES:**

Color & Form Brownish solid
Purity ≥ 95% by TLC
Solubility Soluble in DMF
Absorption/Emission
Extinction
Coefficient

STORAGE AND HANDLING: Store at ≤ -20°C

APPLICATION: *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, complexing of the chelator with copper can be used to selectively cleave RNA or DNA in the presence of hydrogen peroxide.

Ref.: 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) *Proc. Natl. Acad. Sci. USA* **86**, 9702(1989).

TOXICITY: Not established. Not listed by NTP, IARC or OSHA.

FIRST AID:	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.
-------------------	---

Disclaimer: <i>Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.</i>
--