

# ReadiUse™ Universal Kinase Reaction Stopping Buffer

Catalog Number: 31005

Unit Size: 100 mL

## Product Details

---

Storage Conditions	Refrigerated (2-8 °C)
Expiration Date	12 months upon receiving

## Chemical Properties

---

Appearance	Liquid
Molecular Weight	N/A
Soluble In	N/A

## Spectral Properties

---

Excitation Wavelength	N/A
Emission Wavelength	N/A

## Applications

---

ReadiUse™ Universal Kinase Reaction Stopping Buffer is an aqueous solution that effectively chelates divalent metal ions. It is prepared by dissolving Na<sub>2</sub>EDTA•2H<sub>2</sub>O in distilled, deionized water and adjusting the pH to 8.0 using sodium hydroxide. This sterile-filtered solution finds wide applicability in biochemistry and molecular biology research. It is particularly useful for studying cell biology, molecular biology, bioactive small molecules, biochemicals, solutions and reagents, inorganic salts, chelators, and EDTA solutions. Notably, it has been utilized in a study to investigate the role of antigen retrieval in immunostaining ethanol-fixed smear.