

# ReadiUse™ ATP Reference Standard \*10 mM Aqueous Solution\*

Catalog Number: 17198

Unit Size: 5 mL

## Product Details

---

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

## Chemical Properties

---

Appearance	Liquid
Molecular Weight	551.14
Soluble In	Water

## Spectral Properties

---

Excitation Wavelength	N/A
Emission Wavelength	N/A

## Applications

---

The ReadiUse™ ATP Reference Standard \*10 mM Aqueous Solution\* is designed for the accurate quantification of adenosine triphosphate (ATP) in various scientific assays. This standard is formulated with high purity and supplied in an aqueous solution for immediate use in calibrating ATP measurement assays. The ReadiUse™ ATP Reference Standard is applicable for various scientific purposes, including calibration of bioluminescence assays for ATP detection, assessment of microbial contamination, evaluation of cell viability and proliferation, investigation of energy metabolism pathways, and environmental monitoring for microbial load. It is particularly relevant for research and quality control in industries where precise ATP quantification is critical, such as the food and beverage, pharmaceutical, and cosmetic sectors.