

ReadiUse™ APS-5

Catalog Number: 11591

Unit Size: 10 mg

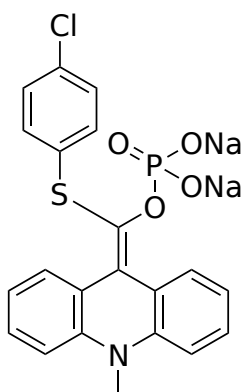
Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	489.82
Soluble In	Water

Chemical Structure



Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

ReadiUse™ APS-5 is a ready-to-use chemiluminescent substrate designed for the sensitive detection of alkaline phosphatase (AP) in ELISA, Western blot, and solution-based assays. Utilizing acridan-based chemistry, ReadiUse™ APS-5 produces a rapid and sustained luminescent signal upon reaction with AP-labeled molecules, enabling highly sensitive and quantitative analysis. The reagent offers excellent performance, with a detection limit below 10^{-19} moles of AP, making it ideal for applications that demand low-level detection and high precision. Its fast signal development significantly shortens assay time and improves throughput, while the extended signal duration provides flexibility in read timing without compromising linearity or accuracy. ReadiUse™ APS-5 exhibits a broad dynamic range with linear calibration across multiple orders of magnitude, and is robust against temperature variations, maintaining consistent performance between 22°C and 35°C. This eliminates the need for strict temperature control during routine assays.