

AMPPD Substrate

Catalog Number: 26111

Unit Size: 1 mg

Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	382.35
Soluble In	Water

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

AMPPD Substrate is a dioxetane-based chemiluminescent substrate used for the detection of alkaline phosphatase (AP) activity in biochemical and immunological assays. Upon enzymatic dephosphorylation, the substrate forms an unstable intermediate that undergoes decomposition, resulting in light emission. This chemiluminescent reaction enables detection of enzyme-labeled targets without the need for external excitation. The emission is glow-type, providing a sustained signal suitable for flexible signal acquisition.

The substrate is commonly used in applications such as Western blotting, ELISA, and other AP-based assays where stable signal output is required. The generated chemiluminescent signal can be detected using imaging systems or film, supporting qualitative and quantitative analysis of biomolecular targets.