

Lumigen APS-5 Substrate

Catalog Number: 26110

Unit Size: 5 mg

Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	489.82
Soluble In	Water

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

Lumigen APS-5 Substrate is a chemiluminescent substrate for detection of alkaline phosphatase (AP) activity in biochemical and immunological assays. Upon enzymatic dephosphorylation by alkaline phosphatase, the substrate undergoes a chemical reaction that produces light, allowing sensitive detection of labeled targets without the need for external excitation.

The substrate is commonly used in applications such as Western blotting, ELISA, and other AP-based assays where high sensitivity and low background are required. Its chemiluminescent signal can be captured using imaging systems or film, supporting both qualitative and quantitative analysis of protein expression and biomolecular interactions.