

CSPD Substrate

Catalog Number: 26112

Unit Size: 5 mg

Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	460.75
Soluble In	Water

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

CSPD Substrate is a chemiluminescent alkaline phosphatase substrate designed for detection of alkaline phosphatase and alkaline phosphatase-labeled molecules in both membrane-based and solution-based assays. Upon enzymatic dephosphorylation, the substrate undergoes a chemical reaction that produces light, enabling sensitive detection without external excitation. Maximum light output is typically reached within minutes, followed by sustained glow-type emission that persists for extended periods.

The substrate is suitable for applications including Western, Southern, and Northern blotting on nitrocellulose membranes, as well as solution-based assays such as immunoassays, DNA probe assays, enzyme assays, and reporter gene assays. It provides low background luminescence and consistent signal output, supporting reliable detection of labeled targets. Chemiluminescent enhancers may be used to increase signal intensity and sensitivity depending on assay format.