

CDP Brite Substrate

Catalog Number: 26113

Unit Size: 1 mg

Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	495.2
Soluble In	Water

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

CDP Brite Substrate is a chemiluminescent alkaline phosphatase substrate used for detection of alkaline phosphatase and alkaline phosphatase-labeled molecules in biochemical and molecular biology assays. Upon enzymatic dephosphorylation, the substrate undergoes a chemical reaction that produces light, enabling detection of labeled targets without external excitation. Maximum light output is typically reached within several minutes, followed by sustained emission that allows flexible signal acquisition.

Chemiluminescent substrates provide a sensitive chemiluminescent alternative to commonly used fluorogenic substrates such as methylumbelliferyl phosphate (MUP) and colorimetric substrates such as p-nitrophenyl phosphate (pNPP). The substrate is suitable for membrane-based applications including Western, Southern, and Northern blotting, as well as solution-based assays such as immunoassays, DNA probe assays, enzyme assays, and reporter gene assays.