

DPX

DPX is a positively charged quencher that is often used with ANTS ([90010](#)) to study membrane fusion or permeability.



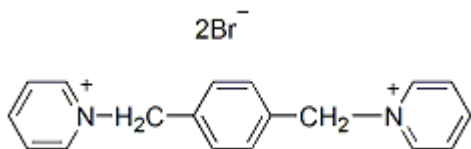
Product attributes

CAS number	14208-10-7
Probe cellular localization	Fluid phase tracer
Cell permeability	Membrane impermeant

Product Description

DPX is a positively charged quencher that is often used with ANTS ([90010](#)) to study membrane fusion or permeability.

- Off-white solid soluble in water
- Store desiccated at 4 °C
- C₁₈H₁₈Br₂N₂
- MW: 422
- [14208-10-7]



This datasheet was generated on June 19, 2026 at 04:45:30 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/dpx/>