



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

## DPX

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



### Product attributes

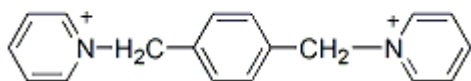
CAS number	14208-10-7
Probe cellular localization	Fluid phase tracer
Cell permeability	Membrane impermeant

## 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
- C<sub>18</sub>H<sub>18</sub>Br<sub>2</sub>N<sub>2</sub>
- MW: 422
- [14208-10-7]

2Br<sup>-</sup>



This datasheet was generated on January 14, 2026 at 05:36:05 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/dpx/>