

## **DPX**

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



Call us: 800-304-5357

## **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

This datasheet was generated on November 29, 2025 at 01:30:04 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/dpx/