## **SDIP/Europium for Membrane Fusion Assay**

stétum

Product attributes

Call us: 800-304-5357

Probe cellular localization	Fluid phase tracer
Cell permeability	Membrane impermeant
Colors	Far-red
Excitation/Emission	250 -320/610 nm

Biotium developed SDIP/Eu<sup>3</sup> as an alternative to DPA/Tb<sup>3</sup> (80104), which is used for vesicle fusion assays.

## **Product Description**

Biotium developed SDIP/Eu³ as an alternative to DPA/Tb³ (80104), which is used for vesicle fusion assays. Similar to DPA/Tb³ assay, one population of vesicles can be loaded with the chelator SDIP while another population of vesicles can be loaded with EuCl₃. Intense red fluorescence is produced when the two types of vesicles fuse as a result of formation of the SDIP/Eu³ complex. The main advantage of SDIP/Eu³ over DPA/TB³ is that the SDIP/Eu³ generates fluorescent emission with brighter intensity and much longer wavelength. Each set of products contain 50 mg SDIP and 25 mg EuCl₃ in two separate vials.

- $\lambda Abs/\lambda Em(for complex) = 250 \sim 320/\sim 610 nm$
- SDIP is a light yellow solid
- EuCl<sup>3</sup> is in a colorless crystal
- Both are readily soluble in water
- Both components are stable at room temperature or 4°C
- Aqueous solutions of SDIP should be protected from light
- MW of SDIP: 449
- MW of EuCl<sup>3</sup>: ~258

This datasheet was generated on November 29, 2025 at 12:58:18 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/sdipeuropium-for-membrane-fusion-assay/">https://biotium.com/product/sdipeuropium-for-membrane-fusion-assay/</a>