

46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

www.biotium.com

PRODUCT AND SAFETY DATA SHEET

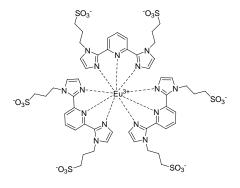
PRODUCT NAME: SDIP/Europium for membrane fusion assay

CATALOG # 80105

COMPONENT Component A: 50 mg SDIP

Component B: 25 mg EuCl₃

MOLECULAR MWt of SDIP: 449 **INFORMATION:** Mwt. of EuCl₃: ~258



PROPERTIES:

Color & Form SDIP is a light yellow solid.

EuCl₃ is in a colorless crystal form.

Solubility Both components are readily soluble in H_2O .

Absorption/Emission $\lambda_{abs} = 250-320 \text{ nm (for complex)}; \lambda_{em} \sim 610 \text{ nm (for complex)}$

STORAGE AND

Both components are stable at room temperature or 4°C. Aqueous solution of SDIP

HANDLING: should be protected from light.

APPLICATION: SDIP/Europium can be used for vesicle fusion assays, similar to the use of DPA/Tb³⁺

(Nature **281**, 690(1979); Biochemistry **19**, 6011(1980); Biochemistry **33**, 5805(1994); J. Biol. Chem. **269**, 14473(1994)). Neither the ligand SDIP nor Eu3⁺ is fluorescent in water. However, when SDIP and Eu³⁺ are combined at 3 to 1 or greater a ratio strong red fluorescence forms due to formation of SDIP/Eu³⁺ complex. High concentrations of phosphate, amino acids, or citrate will interfere with the complex formation and thus should be avoided. We recommend one population of vesicles be loaded with ~0.2mM EuCl₃ and the other population of vesicles be loaded with 1-2 mM SDIP. Including Ca²⁺ and EDTA in the external medium inhibits fluorescent complex formation outside the fused vesicles. Fluorescence is collected at ~610nm, with excitation at 250-320 nm.

TOXICITY: Not established. Not listed by NTP, IARC or OSHA.

FIRST AID: Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with

plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing,

move individual to fresh air and seek medical advice immediately.

Disclaimer: Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.