

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:** DiIC<sub>1</sub>(5) (1,1'-dimethyl-3,3,3',3'-tetramethylindodicarbocyanine, iodide)

CATALOG NUMBER: 70015

**MOLECULAR**  $C_{27}H_{31}IN_2$  **INFORMATION:** Mwt: 510.46

[36536-22-8]

 $H_3C$   $CH_3$   $H_3C$   $CH_3$   $CH_3$ 

**PROPERTIES:** 

Color & FormGreen solidPurity $\geq$  99% by TLCSolubilitySoluble in DMSO $\lambda_{ex} \mid \lambda_{em}$ 646/666nm

**Extinction** 226,000(646 nm; in DMSO)

Coefficient  $(\epsilon)$ 

STORAGE AND HANDLING:

Store desiccated at  $\leq$  4 °C and protect from light, especially in solution.

**APPLICATION:** The carbocyanine dye DiIC<sub>1</sub>(5) has been used to measure mitochondrial membrane

potentials in apoptotic cells. Loss of membrane potential in apoptotic cells is reflected

by a decrease of fluorescence signal in the infrared channel.

Ref: 1) Cytometry 33, 333(1998).

TOXICITY: Unknown

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.