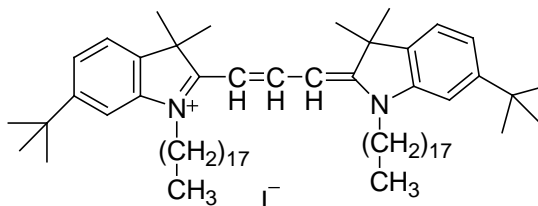


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Neuro-DiI in vegetable oil**CATALOG #:** 60020**MOLECULAR INFORMATION:** C₆₇H₁₁₅N₂
Mwt: 1072**PROPERTIES:**

Color & Form Dark orange oil
Purity ≥ 95% by TLC
Absorption\Emission 549 nm \ 565 nm (in MeOH)
Extinction Coefficient 148,000 cm⁻¹ M⁻¹ (in MeOH)

STORAGE AND HANDLING:Store at room temperature and protect from light. **DO NOT FREEZE.****APPLICATION:**

Neuro-DiI is a proprietary product of Biotium. It is a derivative of DiI and has spectral property similar to that of the later. Neuro-DiI is uniformly dissolved in vegetable oil in a formulation optimized for microinjection. The green fluorescent Neuro-DiO (60019) and red fluorescent DiI in oil can be used in combination for two color imaging.

It is recommended to heat-vortex **60020** at least 10 min at 55C each time prior to use.

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.