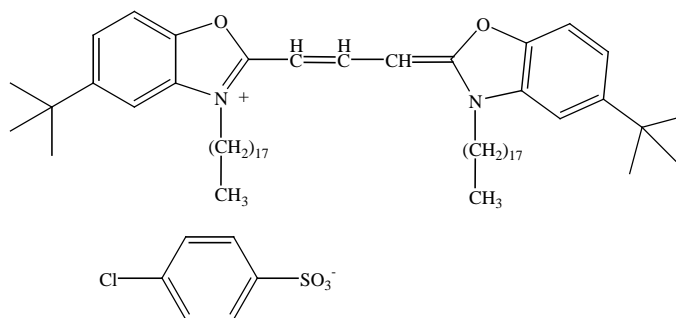


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Neuro-DiO in vegetable oil**CATALOG #:** 60019**MOLECULAR INFORMATION:** C₆₇H₁₀₅ClN₂O₅S
Mwt: 1086**PROPERTIES:**

Color & Form Yellow orange oil
Dye Concentration 0.25% of the dye by weight in the oil (density of oil ~0.8mg/uL)
Purity ≥ 95% by TLC
Absorption\ Emission 484 nm \ 501 nm (in MeOH)
Extinction Coefficient 150,000 cm⁻¹ M⁻¹ (in MeOH)

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

Neuro-DiO is a proprietary product of Biotium. It is a derivative of DiO and has spectral property similar to that of the later. However, unlike DiO, Neuro-DiO is structurally more similar to DiI, which is known to be a much better lipophilic neuronal tracer than regular DiO. Like DiI in oil, Neuro-DiO is uniformly dissolved in vegetable oil in a formulation optimized for microinjection. The green fluorescent Neuro-DiO and red fluorescent DiI in oil can be used in combination for two color imaging.

It is recommended to heat-vortex **60019** 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.
