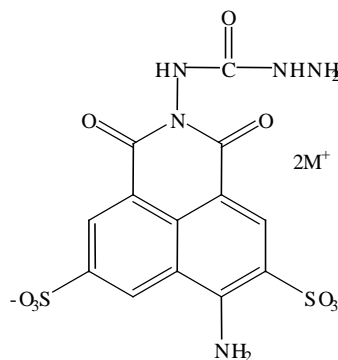


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** Lucifer Yellow CH lithium salt, 100mM in water**CATALOG #** 80026**MOLECULAR INFORMATION:** C<sub>14</sub>H<sub>10</sub>Li<sub>2</sub>N<sub>5</sub>O<sub>9</sub>  
S<sub>2</sub>  
MWt: 45780015 M = Li<sup>+</sup>  
80016 M = K<sup>+</sup>**PROPERTIES:****Color & Form** Light orange solution  
**Purity** ≥ 95% by HPLC  
**Absorption/Emission** 428/536 nm**STORAGE AND HANDLING:** Stored at room temperature and protect from light.**APPLICATION:** Lucifer Yellow CH lithium salt (LY CH lithium salt) is a widely used polar molecular tracer for studying neuronal morphology. The fluorescent molecule contains a carbohydrazide that allow the molecule to be fixable with an aldehyde fixation agent. The solution of Lucifer Yellow CH is commonly used for microinjection.Ref.: 1) *Nature* **292**, 17(1981); 2) *Cell* **14**, 741(1978).**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.