

## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** CF640R-UTP, Lyophilized powder

**CATALOG #** 40032

**PROPERTIES:**

**Color & Form** Blue solids  
**Solubility** Soluble in water

**STORAGE AND HANDLING:** Store desiccated at -20°C, protect from light.

**APPLICATION:** CF640R-UTP is a fluorescent labeled UTP. It can be enzymatically incorporated into probes for use in many biological applications. Far-red fluorescent CF640R is spectrally similar to Alexa Fluor 647 but is much more photostable

**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.*