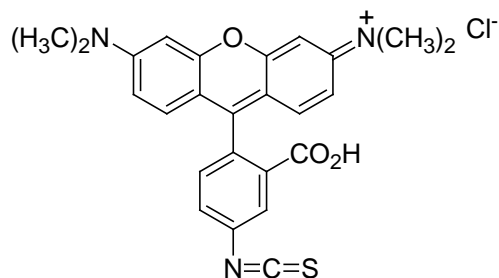


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** Tetramethylrhodamine-5-isothiocyanate isomer G (5-TRITC, G isomer)**CATALOG #** 90018**MOLECULAR INFORMATION:** C<sub>25</sub>H<sub>22</sub>ClN<sub>3</sub>O<sub>3</sub>S  
Mwt: 479.98**PROPERTIES:**

**Color & Form** Dark red solid  
**Purity** ≥ 90% by TLC  
**Solubility** Soluble in DMF  
**Absorption/Emission** 544nm/572nm (MeOH)  
**Extinction Coefficient** >75,000

**STORAGE AND HANDLING:** Store desiccated at -20 °C. Protect from light, especially when in solution.**APPLICATION:** 5-TRITC is spectrally similar to TAMRA. The isothiocyanate reactive group reacts with primary or secondary amines to form a thiourea bond linkage.**RELATED PRODUCTS** *Tetramethylrhodamine-6-isothiocyanate isomer R (6-TRITC, R isomer) (#90019);  
5-Carboxytetramethylrhodamine-PEO<sub>8</sub>-propionate succinimidyl ester(#90087);  
5-Carboxytetramethylrhodamine-PEO<sub>12</sub>-propionate succinimidyl ester(#90088)***TOXICITY:** Not established. Not listed by NTP, IARC or OSHA.

<b>FIRST AID:</b>	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.
-------------------	---

<b>Disclaimer:</b> <i>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.</i>
--