

46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352 www.biotium.com

## **PRODUCT AND SAFETY DATA SHEET**

**PRODUCT NAME:** 

TMP-PEG3-amine, TFA salt

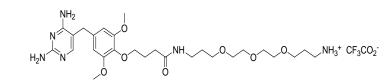
CATALOG #

**MOLECULAR** 

**INFORMATION:** 

C<sub>29</sub>H<sub>45</sub>F<sub>3</sub>N<sub>6</sub>O<sub>9</sub> MWt: 678.71

91056



| PROPERTIES:<br>Color & Form<br>Solubility<br>STORAGE AND<br>HANDLING: | White solid<br>Soluble in DMF or DMSO<br>Store at -20°C   |
|---|---|
| APPLICATION:  | The trimethoprim-PEG3-amine (TMP-PEG3-amine) can be conjugated to various probes to make trimethoprim-based chemical tag (TMP-tag). The TMP-tag is based on the high affinity ( $K_D \sim 1nM$ ) binding of TMP to the <i>E.coli</i> dihydrofolate reductase (eDHFR). The target protein, fused to eDHFR via standard molecular cloning technology, is labeled by the TMP-probe in a non-covalent manner. The low affinity of TMP with mammalian DHFR ( $K_D > 1\mu M$ ) ensures minimal background. The TMP-tag has been demonstrated to enable protein visualization and manipulation both inside and on surface of mammalian cells.[1-4] |
|   | Ref.: 1) Nat. Methods <b>2</b> , 255(2005); 2) ChemBioChem <b>8</b> , 767(2007); 3) Nat. Methods <b>7</b> , 717(2010); 4) J Am Chem Soc. <b>133</b> , 10006(2011)   |
| TOXICITY:   | Not established. Not listed by NTP, IARC or OSHA.   |
| 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.*