

TMP-PEG3-Amine, TFA salt

The trimethoprim-PEG3-amine (TMP-PEG3-amine) can be conjugated to various probes to make trimethoprim-based chemical tag (TMP-tag).



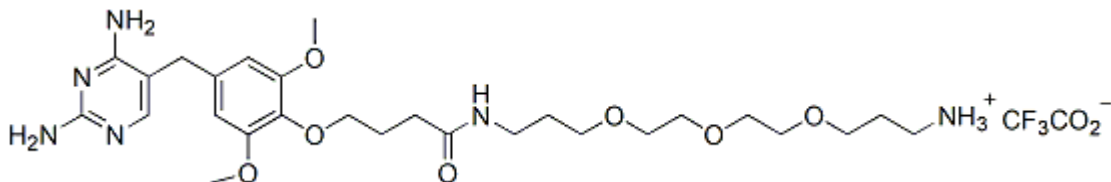
Product attributes

Product Description

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 ($K_d \sim 1$ nM) binding of TMP to the *E. coli* 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 ($K_d > 1$ uM) ensures minimal background. The TMP-tag has been demonstrated to enable protein visualization and manipulation of both intracellular and surface targets in mammalian cells (1-4).

The TMP-PEG-amine can be used in combination with our [Mix-n-Stain™ Small Ligand Labeling Kits](#) to generate fluorescent chemical tags without purification for cell surface or intracellular staining.

- White solid soluble in DMF or DMSO
- MW: 678.71



References

1. Miller, L.M et al (2005) Nat. Methods 2:255-257.
2. Calloway, NT. et al (2007) ChemBioChem 8:767-774.
3. Wombacher, R. et al (2010) Nat. Methods 7:717-719
4. Long, MJ. et al (2011) J Am Chem Soc. 133:10006–10009

This datasheet was generated on January 8, 2026 at 04:05:11 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/tmp-peg3-amine-tfa-salt/>