



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

## Biotin-PEO2-PPO2-amine, TFA salt

Biotin-PEO<sub>2</sub>-PPO<sub>2</sub>-amine trifluoroacetic acid (TFA) salt is a water soluble biotin derivative with an amine group.

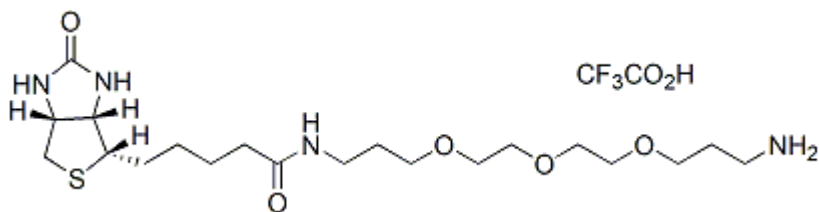


### Product attributes

## Product Description

Biotin-PEO<sub>2</sub>-PPO<sub>2</sub>-amine trifluoroacetic acid (TFA) salt is a water soluble biotin derivative with an amine group.

- White solid soluble in water or DMSO
- Store at 4 °C
- C<sub>22</sub>H<sub>39</sub>F<sub>3</sub>N<sub>4</sub>O<sub>7</sub>S
- MW: 560.63



This datasheet was generated on December 22, 2025 at 12:18:51 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/biotin-peo2-ppo2-amine-trifluoroacetic-acid-salt/>