Biotin-PEO2-PPO2-amine, TFA salt

Biotin-PEO₂-PPO₂-amine trifluoroacetic acid (TFA) salt is a water soluble biotin derivative with an amine group.



Product attributes

Product Description

Biotin-PEO₂-PPO₂-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
- C₂₂H₃₉F₃N₄O₇S
- MW: 560.63

This datasheet was generated on November 3, 2025 at 10:57:44 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/biotin-peo2-ppo2-amine-trifluoroacetic-acid-salt/