

## Biotin NTA

Biotin NTA, also known as BNTA (Biotin nitrilotriacetic acid, tripotassium salt) is used to detect histidine-tagged proteins (as little as 0.1 pmol) immobilized on nitrocellulose membranes.

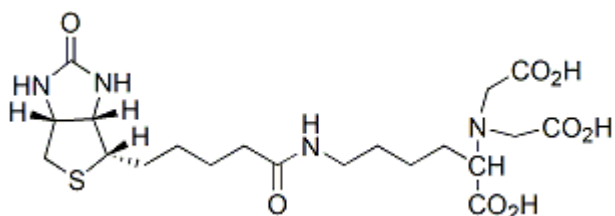


### Product attributes

## Product Description

Biotin NTA, also known as BNTA (Biotin nitrilotriacetic acid, tripotassium salt) is used to detect histidine-tagged proteins (as little as 0.1 pmol) immobilized on nitrocellulose membranes. Biotin NTA can be loaded with Ni by combining it with a 3 fold molar excess of  $\text{NiSO}_4$  at pH7.9 (1). Biotin NTA can be removed from the histidine-tagged protein at pH 4.8, allowing the blot be reanalyzed with another probe.

- White solid soluble in water
- Store at 4°C
- $\text{C}_{20}\text{H}_{36}\text{N}_4\text{O}_8\text{S}$
- MW: 488.56



## References

1. Anal Biochem 236, 101 (1996).
2. Anal Biochem 229, 119 (1995).

This datasheet was generated on December 22, 2025 at 11:57:39 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/biotin-nta-biotin-nitrilotriacetic-acid-tripotassium-salt-also-called-bnta/>