Call us: 800-304-5357 Email: btinfo@biotium.com

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 NiSO4 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
- C₂₀H₃₆N₄O₈S
- MW: 488.56

References

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

This datasheet was generated on November 5, 2025 at 05:44:27 PM. 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/