

Biotin-X NTA

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

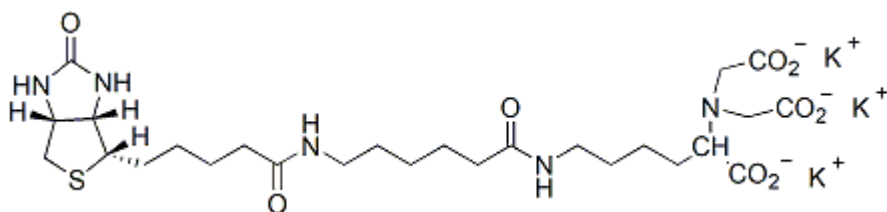


Product attributes

Product Description

Biotin-X NTA (Biotin-X nitrilotriacetic acid, tripotassium salt) is used to detect histidine-tagged proteins (as little as 0.1 pmol) immobilized on nitrocellulose membranes. Biotin-X NTA can be loaded with Ni by combining it with a 3 fold molar excess of NiSO₄ at pH7.9(1). Biotin-X NTA can be removed from the histidine-tagged protein at pH 4.8, allowing the blot to be reanalyzed with another probe.

- White solid soluble in water
- Store at 4°C
- C₂₆H₄₀K₃N₅O₉S
- MW: 715.98



References

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

This datasheet was generated on June 3, 2026 at 07:14:38 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/biotin-x-nta-or-biotin-x-nitrilotriacetic-acid-tripotassium-salt/>