

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.

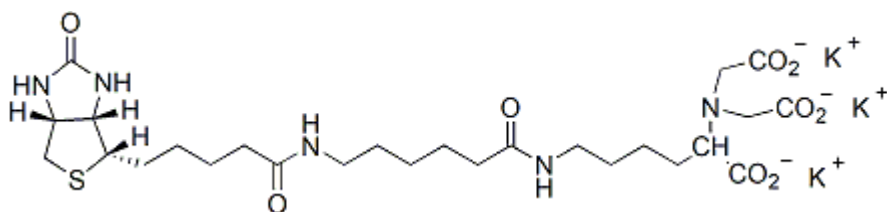


Product attributes

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 NiSO_4 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
- $\text{C}_{26}\text{H}_{40}\text{K}_3\text{N}_5\text{O}_9\text{S}$
- MW: 715.98



References

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

This datasheet was generated on December 19, 2025 at 06:31:33 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/>