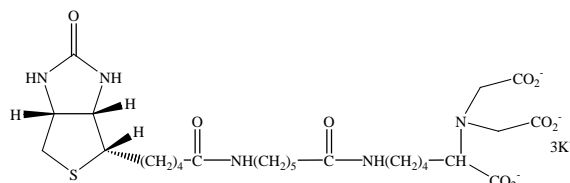


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: **Biotin-X NTA** (biotin-X nitrilotriacetic acid, tripotassium salt)**CATALOG #:** **90056****MOLECULAR INFORMATION:** C₂₆H₄₀K₃N₅O₉S
MW: 715.98**PROPERTIES:**

Color & Form White solid
Purity ≥ 95% by TLC
Solubility Soluble in water

STORAGE AND HANDLING: Store at 4°C**APPLICATION:** Biotin-X NTA is used to detect histidine-tagged proteins (for as little as less than 0.1 pmol) immobilized on nitrocellulose membranes. Biotin-X NTA can be removed from the histidine-tagged protein at pH 4.8, allowing the blot to be reanalyzed with another probe.Ref.: 1) *Anal. Biochem.* **236**, 101 (1996); 2) *Anal. Biochem.* **229**, 119 (1995).**TOXICITY:** Unknown.

FIRST AID:	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.
-------------------	---

Disclaimer: <i>Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.</i>
--