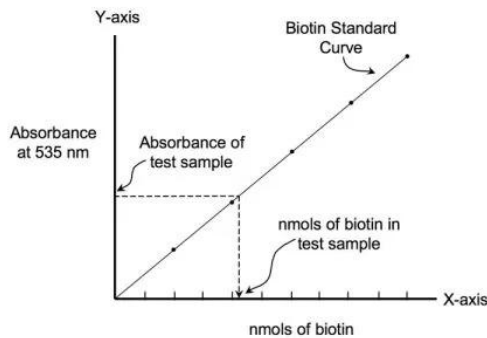




QUANTTAG BIOTIN QUANTITATION KIT

SKU: BDK-2000



DESCRIPTION

The QuantTag Biotin Kit (BDK-2000) is designed to determine the amount of free biotin in solution or the number of biotin molecules attached to nucleic acids, proteins, or other macromolecules. This kit can be used to accurately determine the labeling efficiency of your process to label molecules with biotin.

SPECIFICATIONS

Unit Size 1 Kit

Applications Immunohistochemistry / Immunocytochemistry, Immunofluorescence, In situ hybridization, Blotting Applications, Elispot, ELISAs, Flow Cytometry/Cell Separation, Glycobiology

Safety Title WARNING: NA -www.P65Warnings.ca.gov

For research use only. Not intended for therapeutic or diagnostic use in animals or humans.



KIT COMPONENTS

- Reagent 1 (87.5 ml)*
- Reagent 2 (87.5 ml)*
- Reagent 3 (8.75 ml)*
- Biotin Standard Solution (1 ml; 0.5 mM)

The kit contains enough material to perform 25 assays using the 1 ml protocol (including the creation of a standard curve for each) or 250 assays using the 0.1 ml protocol.

*Note: Avoid contact with skin and eyes. If contact occurs, flush immediately with water.

TECHNICAL INFORMATION

The QuantTag Biotin Kit (BDK-2000) is designed to determine the amount of free biotin in solution or the number of biotins attached to nucleic acids, proteins, or other macromolecules. This kit can be used to accurately determine the labeling efficiency of your process.

Unlike conventional biotin quantitation methods like the HABA assay, no predigestion of protein or nucleic acids is required, which saves time and increases accuracy. The QuantTag Biotin Kit, being more sensitive than the HABA assay, is able to detect less than 1 nmol of biotin.

QuantTag Kit reagents chemically react with free or bound biotin producing a colored product that can be quantified using a spectrophotometer. The absorbance is measured in the visible spectrum allowing the use of plastic cuvettes or microtitre plates.

The QuantTag Biotin Kit is quick and easy to use and the assay can be completed in 30 minutes. The kit contains sufficient material to perform from 25 to 250 tests depending on assay size. A biotin standard is included.

CITATIONS



Powered by Bioz © 2023 See more details on Bioz

DOCUMENTS

- [User Guide](#)

For research use only. Not intended for therapeutic or diagnostic use in animals or humans.



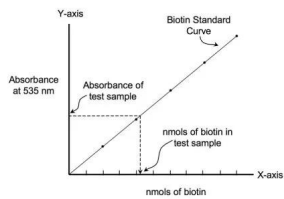
www.vectorlabs.com

Email: customerservice@vectorlabs.com

Telephone: [\(650\) 697-3600](tel:(650)697-3600)

- [Safety Data Sheet](#)
- [Download CoA](#)
- [Datasheet](#)

GALLERY IMAGES



For research use only. Not intended for therapeutic or diagnostic use in animals or humans.

[QuantTag Biotin Quantitation Kit](#)

<https://vectorlabs.com/products/quant-tag-biotin-quantitation-kit/>