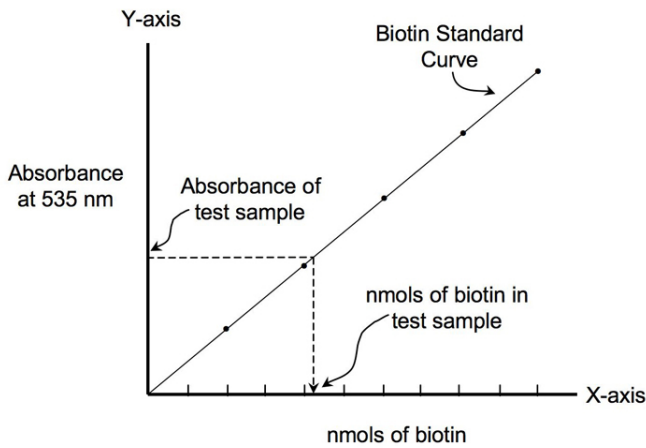




QuantTag Biotin Quantitation Kit

BDK-2000

Product Images



Short 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.

Additional Information

Unit Size	1 kit
Applications	Immunohistochemistry / Immunocytochemistry, Immunofluorescence, In situ hybridization, Blotting Applications, Elispot, ELISAs, Flow Cytometry/Cell Separation, Glycobiology
Safety Title	NA

