



CLICK-&-GO® PROTEIN ENRICHMENT KIT FOR CAPTURE OF ALKYNE-MODIFIED PROTEINS

SKU: CCT-1039



DESCRIPTION

Click-&-Go™ Protein Enrichment Kit is an efficient tool for capture of alkyne-modified proteins on an azide-agarose resin. The alkyne modification can occur via metabolic feeding, enzymatic addition, or chemical modification. The alkyne-modified proteins, or their post-translationally modified forms, are captured from complex protein extracts on the azide-agarose resin supplied. Once covalently attached to the resin via copper catalyzed click chemistry, beads can be washed with highest stringency virtually eliminating any non-specifically bound proteins to yield a highly enriched population of nascent molecules. Upon protease digestion, this yields a highly pure peptide pool that is ideal for mass spectrometry (e.g., LC MS/MS) based analysis.

The biotin-free enrichment improves signal to noise by eliminating nonspecific binding and increasing selectivity, thus improving detection of low abundant proteins. This approach is fully compatible with widely used MS techniques including iTRAC and ICAT.

The biotin-free, covalent capture approach is superior to biotin or lectin based enrichment approaches, and ideal for proteomics, biomarker discovery, posttranslational modification (PTM) analysis, and more.

The Click-&-Go™ Protein Enrichment Kit can be used as direct replacement of Click-iT® Protein

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



Enrichment Kit

SPECIFICATIONS

Enrichment Target	Alkyne-modified proteins
Format	Spin Column
Isolation Technology	Biotin-free covalent capture
Unit Size	1 Kit
Storage Instructions	4C
Purchase Note	<p>Add some note</p>
Shipping Instructions	Ambient temperature

SCHEMATICWORKFLOW

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



www.vectorlabs.com

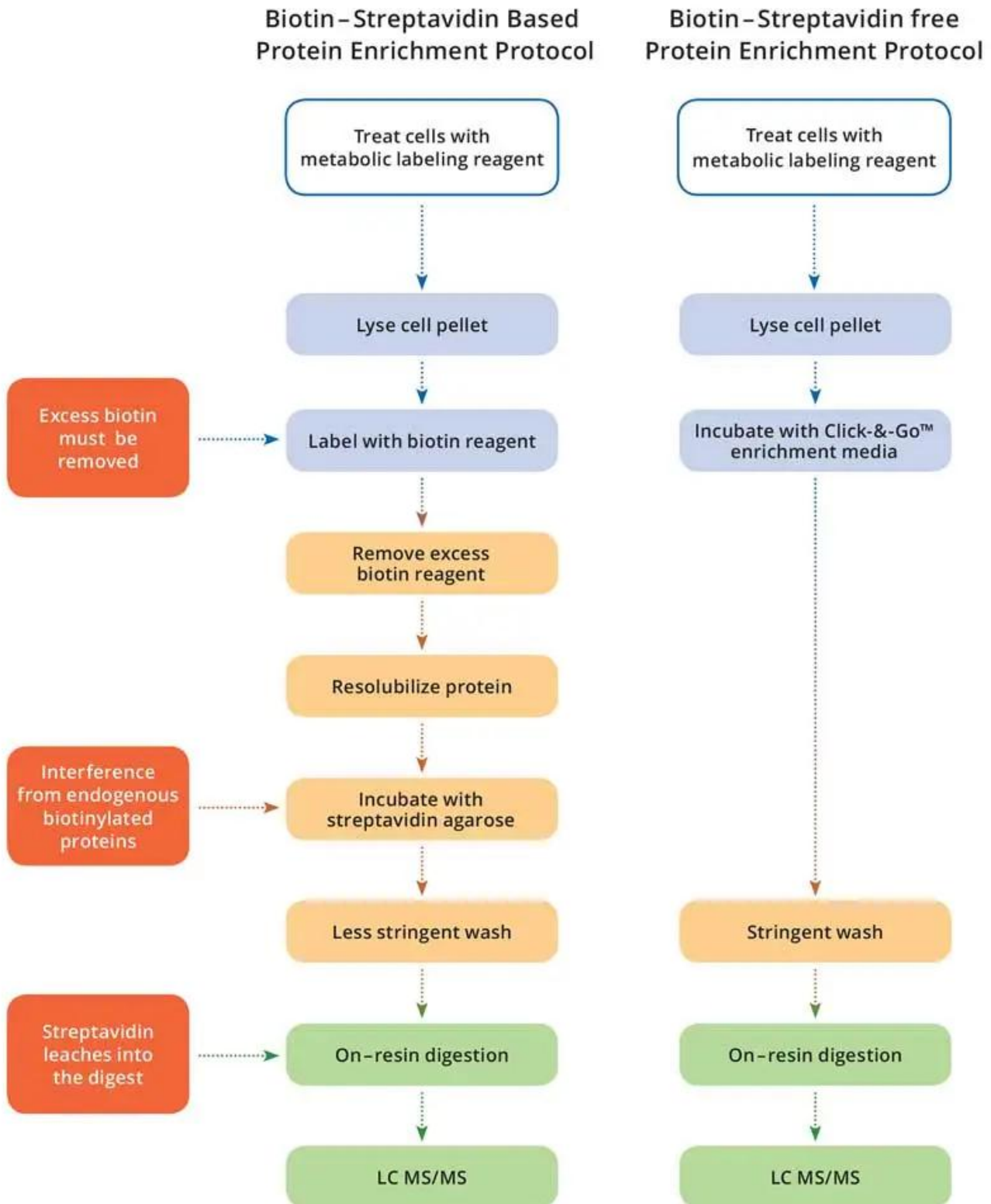
Email: customerservice@vectorlabs.com

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

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

[Click-&-Go® Protein Enrichment Kit for capture of alkyne-modified proteins](https://vectorlabs.com/products/click-go-protein-enrichment-kit-2/)

<https://vectorlabs.com/products/click-go-protein-enrichment-kit-2/>



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



DOCUMENTS

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

GALLERY IMAGES



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