



# **CLICK-&-GO® PROTEIN ENRICHMENT KIT FOR CAPTURE OF AZIDE-MODIFIED PROTEINS**

**SKU:** CCT-1033



---

## **DESCRIPTION**

Click-&-Go® Protein Enrichment Kit is an efficient tool for capture of azide-modified proteins on an alkyne-agarose resin. The azido-modified proteins, or their post-translationally modified forms, are captured from complex protein extracts on the alkyne-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.

## **SPECIFICATIONS**

<b>Enrichment Target</b>	Azide-modified proteins
<b>Format</b>	Spin Column
<b>Isolation Technology</b>	Biotin-free covalent capture
<b>Unit Size</b>	1 Kit
<b>Storage Instructions</b>	4C

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



[www.vectorlabs.com](http://www.vectorlabs.com)

Email: [customerservice@vectorlabs.com](mailto:customerservice@vectorlabs.com)

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

**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.**

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

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



[www.vectorlabs.com](http://www.vectorlabs.com)

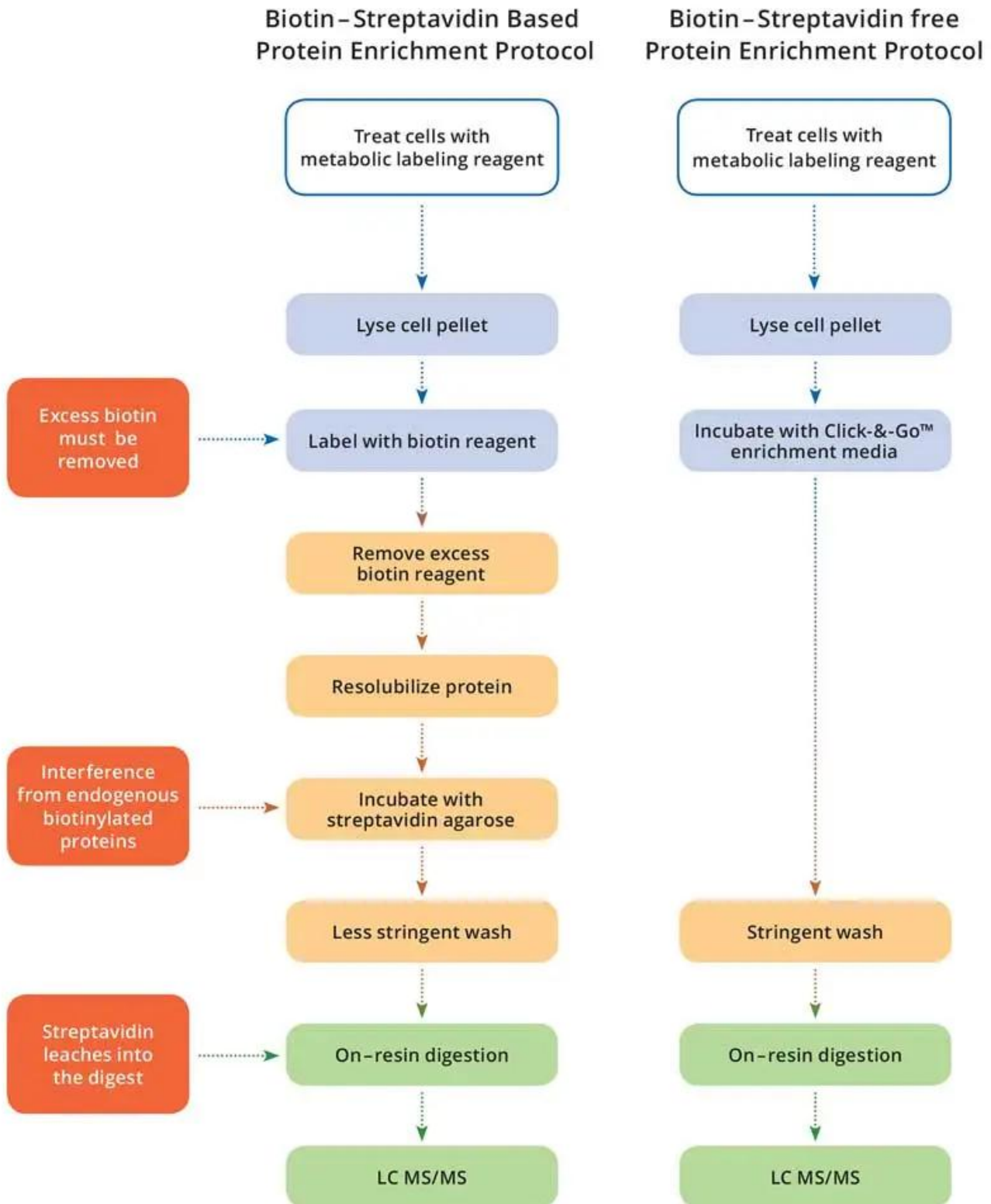
Email: [customerservice@vectorlabs.com](mailto:customerservice@vectorlabs.com)

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 azide-modified proteins](https://vectorlabs.com/products/click-go-protein-enrichment-kit/)

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



**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.**