



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

SKU: CCT-1235



DESCRIPTION

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

SPECIFICATIONS

Enrichment Target	Alkyne-modified proteins
Format	Spin Column
Isolation Technology	Biotin-free covalent capture
Unit Size	1 Kit

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)

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.

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

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



www.vectorlabs.com

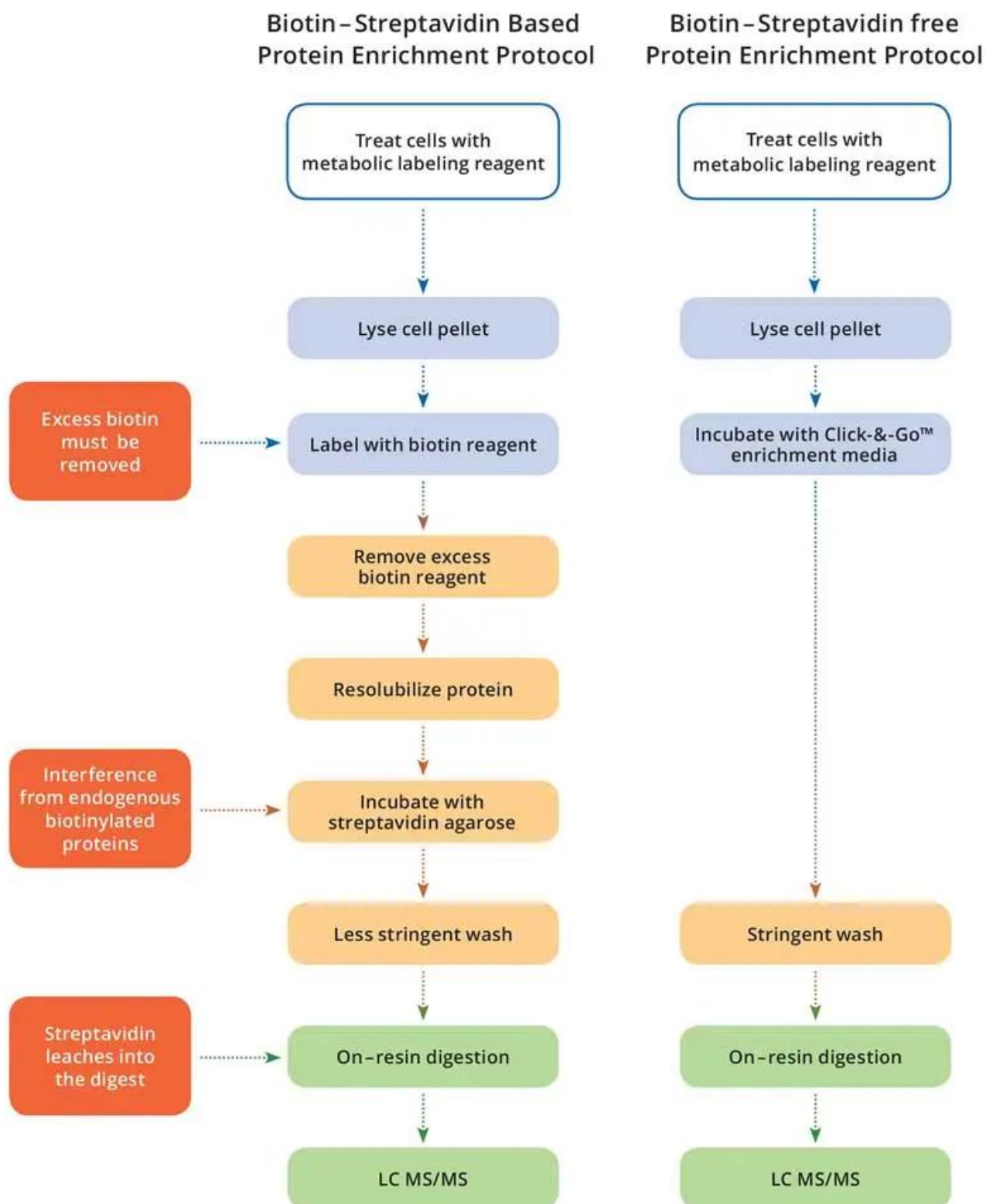
Email: customerservice@vectorlabs.com

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® Plus Protein Enrichment Kit for capture of alkyne-modified proteins](https://vectorlabs.com/products/click-go-plus-protein-enrichment-kit/)

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



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)

DOCUMENTS

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

GALLERY IMAGES



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

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

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