



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

SKU: CCT-1441



DESCRIPTION

Click-&-Go® Protein Capture Kit provides all the necessary reagents to perform capture of alkyne-modified proteins on high capacity streptavidin agarose resin. The kit includes specially formulated components to perform copper-catalyzed click reaction and subsequent capture on high capacity streptavidin agarose. Sufficient material is supplied for 25 enrichments based on the provided protocol. The kit provides alkyne labeled BSA as a positive control.

SPECIFICATIONS

Enrichment Target	Alkyne-modified proteins
Isolation Technology	Biotin-streptavidin based capture
Unit Size	1 Kit
Storage Instructions	4C
Number of Enrichments	25
Purchase Note	Add some note

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)

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 alkyne-modified proteins](https://vectorlabs.com/products/click-go-protein-enrichment-kit-4/)

<https://vectorlabs.com/products/click-go-protein-enrichment-kit-4/>



www.vectorlabs.com

Email: customerservice@vectorlabs.com

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

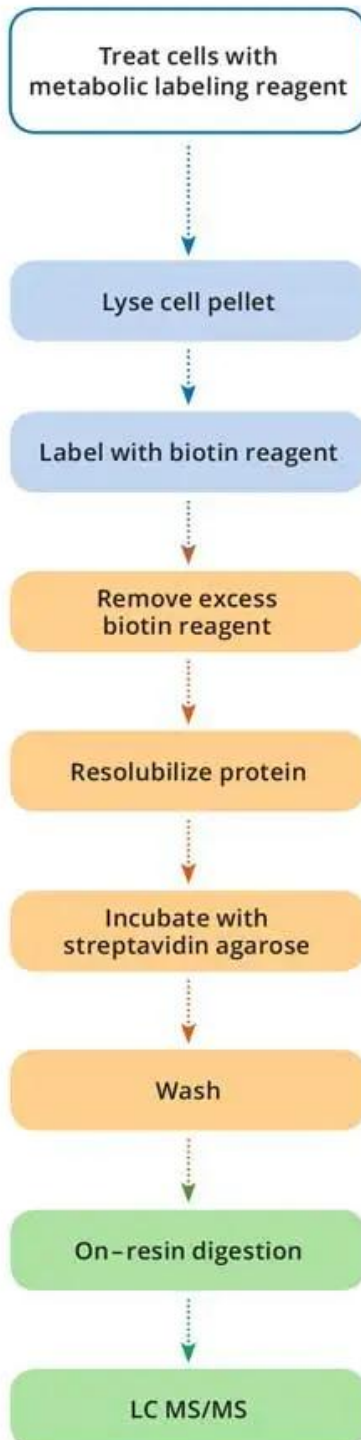
For research use only. Not intended for therapeutic or diagnostic use in animals or humans.

[Click-&-Go® Protein Enrichment Kit for capture of alkyne-modified proteins](https://vectorlabs.com/products/click-go-protein-enrichment-kit-4/)

<https://vectorlabs.com/products/click-go-protein-enrichment-kit-4/>



Biotin–Streptavidin Based Protein Enrichment Protocol



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® Protein Enrichment Kit for capture of alkyne-modified proteins](https://vectorlabs.com/products/click-go-protein-enrichment-kit-4/)

<https://vectorlabs.com/products/click-go-protein-enrichment-kit-4/>