

CLICK-&-GO® PROTEIN ENRICHMENT KIT

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.

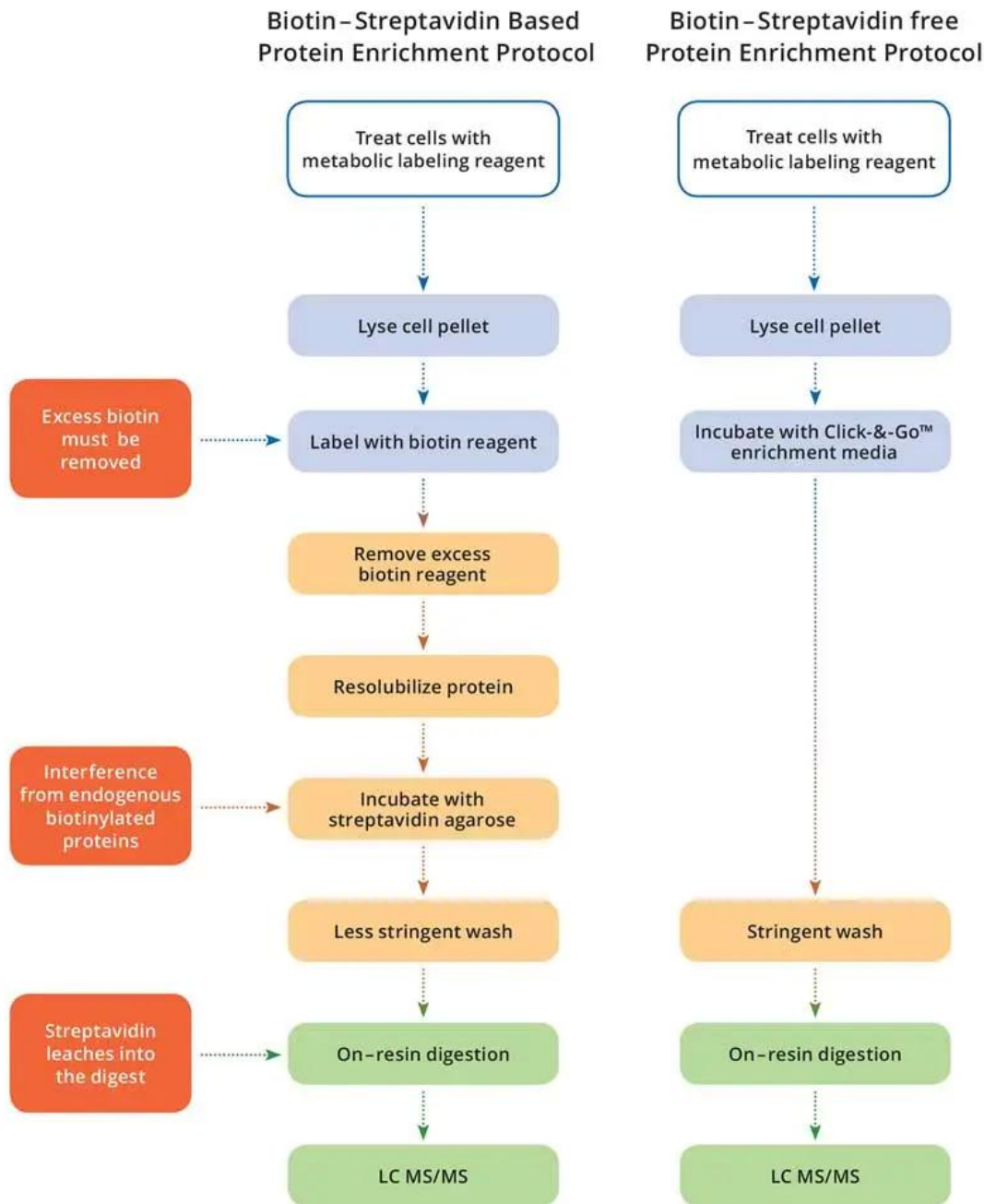
The biotin-free enrichment improves signal to noise by eliminating nonspecific binding and increasing selectivity, thus improving detection of low abundant proteins. This approach is fully compatible with widely used MS techniques including iTRAC and ICAT.

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 Enrichment Kit

Schematic Workflow

For research use only. Not intended for animal or human therapeutic or diagnostic use.



Specifications

Unit Size	1 kit
Enrichment target	Alkyne-modified proteins
Format	Spin Column

For research use only. Not intended for animal or human therapeutic or diagnostic use.

Isolation technology
Storage Conditions
Shipping Conditions

Biotin-free covalent capture
4C
Ambient temperature

For research use only. Not intended for animal or human therapeutic or diagnostic use.