

CLICK-&-GO® PROTEIN ENRICHMENT KIT

SKU: CCT-1441



Description

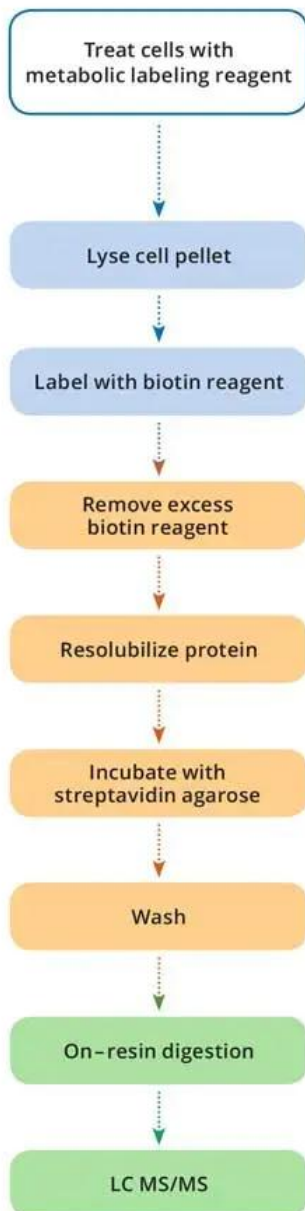
Click-&-Go® Protein Capture Kit provides all the necessary reagents to perform capture of alkyne-modified proteins on high capacity streptavidin agarose resin. The kit includes specially formulated components to perform copper-catalyzed click reaction and subsequent capture on high capacity streptavidin 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
Number of enrichments	25
Isolation technology	Biotin-streptavidin based capture
Storage Conditions	4C
Shipping Conditions	Ambient temperature

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

Biotin-Streptavidin Based Protein Enrichment Protocol



Schematic Workflow

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