

5' Feature Barcode Antibody Conjugation Kit - Lightning-Link® - Oligos 1-10 ab270703

3 Images

Overview

Product name	5' Feature Barcode Antibody Conjugation Kit - Lightning-Link® - Oligos 1-10
Product overview	<p>5' Feature Barcode Antibody Conjugation Kit - Lightning-Link® - Oligos 1-10 (ab270703) enables simple and rapid conjugation of oligonucleotide barcodes to antibodies to enable cell surface protein analysis in Multiomic Cytometry on 10x Genomics Chromium Controller. This kit is compatible with Chromium Single Cell Immune Profiling.</p> <p>The kit includes 10 distinct oligonucleotide barcodes ready to conjugate to 10 µg of antibody each. The protocol involves only a few steps, requires minimal hands-on time and doesn't require any expertise in antibody conjugation or oligonucleotide sourcing.</p> <p>Please find antibody buffer considerations and compatibility in the protocol booklet. For incompatible buffers and low antibody concentrations, please see our antibody purification and concentration kits.</p> <p>For conjugating antibodies with additional distinct oligonucleotide barcodes, please see 5' Feature Barcode Antibody Conjugation Kit - Lightning-Link® - Oligos 11-20 (ab270705) and 5' Feature Barcode Antibody Conjugation Kit - Lightning-Link® - Oligos 21-30 (ab270709). 100 µg conjugation kits are available on-demand for each of these 30 oligonucleotide barcodes. For conjugating more than 30 antibodies with distinct oligonucleotide barcodes, please contact us to discuss your requirements.</p>
Notes	<p>For oligonucleotide barcode sequence information, please see appendix in the protocol booklet.</p> <p>Storing and handling conjugation kits</p> <p>Lyophilized Lightning-Link® components are hygroscopic.</p> <p>Kits are intentionally shipped at ambient temperature with silica gel to avoid exposure to moisture. Upon receipt, store the kit frozen and protect from moisture. Before opening the outer container, allow the lyophilized components to reach room temperature to minimize condensation.</p>

Properties

Storage instructions	Store at -20°C. Please refer to protocols.
-----------------------------	--

Components	1 kit
ab274246 - Conjugate Resuspension Buffer	1 x 1500µl
Modifier reagent	1 x 200µl
ab274247 - Oligo 10X5001	1 x 10µg
ab274248 - Oligo 10X5002	1 x 10µg
ab274249 - Oligo 10X5003	1 x 10µg
ab274250 - Oligo 10X5004	1 x 10µg
ab274251 - Oligo 10X5005	1 x 10µg
ab274252 - Oligo 10X5006	1 x 10µg
ab274253 - Oligo 10X5007	1 x 10µg
ab274254 - Oligo 10X5008	1 x 10µg
ab274255 - Oligo 10X5009	1 x 10µg
ab274256 - Oligo 10X5010	1 x 10µg
ab274244 - Quencher N	1 x 200µl
ab274245 - Reagent C	1 x 500µg
ab274243 - Reagent S	1 x 1250µl

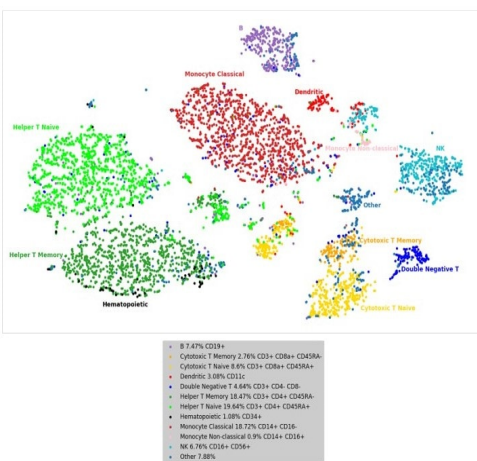
Images



Compatible
Products

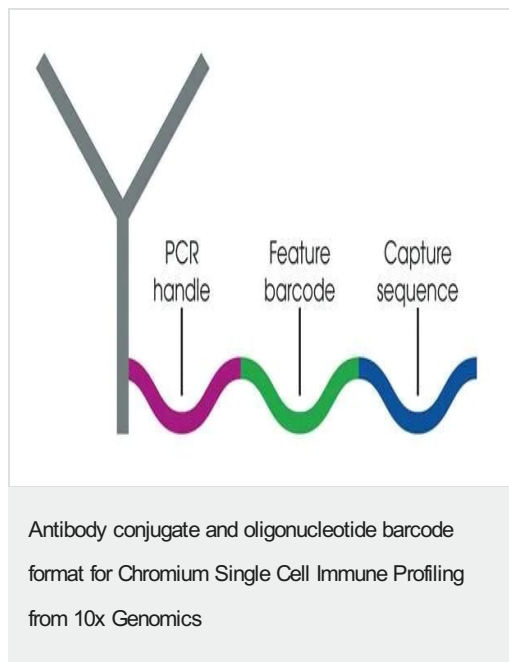
5' Feature Barcode Antibody Conjugation Kit -
Lightning-Link® is part of the 10x Genomics
Compatible Products

5' Feature Barcode Antibody Conjugation Kit - Lightning-Link® is compatible with Chromium Single Cell Immune Profiling from 10x Genomics.



tSNE projections of ~5,000 peripheral blood mononuclear cells (PBMCs) output by Cell Ranger PBMCs were grouped by cell surface protein expression profiles from 10 antibody conjugates for cell surface protein markers generated with 5' Feature Barcode Antibody Conjugation Kit - Lightning-Link®. The additional information from antibody UMI counts better can enrich phenotype information of many canonical immune cell types compared to gene expression data alone - naive and memory T cells, CD4+ and CD8+ T cells (populations denoted with asterisk*), for example.

Cell surface protein analysis using 10 antibody conjugates generated with 5' Feature Barcode Antibody Conjugation Kit - Lightning-Link® on 10x Chromium Controller



5' Feature Barcode Antibody Conjugation Kit - Lightning-Link® includes oligonucleotide barcodes compatible with Chromium Single Cell Immune Profiling from 10x Genomics, which are in the format:

5' CGGAGATGTGTATAAGAGACAGNNNNNNNN
 NNNXXXXXXXXXXXXXXXXNNNNNNNNNCCCAT
 ATAAGA*A*A 3'.

X represents the nucleotides of the unique barcode.

N represents a randomly selected A, C, G, or T.

The symbol * indicates a phosphorothioated bond, this modification renders the internucleotide linkage resistant to nuclease degradation.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors