

TotalSeq™ Universal Cocktails

Instructions for Use

The protocol below is intended for customers who are using TotalSeq™ Universal Cocktails.

Please read the entire protocol below and the appropriate 10x Genomics user guide before starting your experiments. The recommended 10x Genomics user guide for each TotalSeq™ Universal Cocktail is listed in the corresponding TotalSeq™-A, B, or C protocol. BioLegend protocols can be found at: [biolegend.com/en-us/technical-protocols#proteogenomictotalseqprotocols](https://www.biolegend.com/en-us/technical-protocols#proteogenomictotalseqprotocols)

Table 1: Reconstitution and Staining Volumes

TotalSeq™ Format	Reconstitution Volume (μL)	No. of Cells	Volume of FcR Blocked Cells (μL)	Total Staining Volume (μL)
A, B, C	27.5	5x10 ⁵	25	50

Custom cocktails may require different reconstitution instructions. Please reach out to your local BioLegend representative or distributor if you have questions regarding custom cocktails.

Lyophilized Panel Reconstitution and Staining

- Equilibrate the lyophilized panel vial(s) to room temperature for 5 minutes.
- Place lyophilized panel vial in an empty Eppendorf tube, spin down at 10,000 x *g* for 30 seconds at room temperature.
- Rehydrate lyophilized panel by adding 27.5 μL of Cell Staining Buffer (BioLegend Cat. No. **420201**).
Replace the cap and vortex for 10 seconds.
Note: Excess volume added to aid in removal of potential protein aggregates.
- Incubate at room temperature for 5 minutes.
- Vortex again and spin down at 10,000 x *g* for 30 seconds at room temperature.
- Transfer the entire volume (27.5 μL) of reconstituted cocktail to a low protein binding Eppendorf tube (Fisher Cat. No. 022431081 or similar tube).
- Centrifuge at 14,000 x *g* for 10 minutes at 4°C.
- While cocktail is being centrifuged, resuspend and block cells. Perform staining in 12 x 75 mm tubes.
 - For human cells add 2.5 μL of Human TruStain FcX™ Fc blocking reagent (BioLegend Cat. No. **422301**) to 5 x 10⁵ cells in 22.5 μL Cell Staining Buffer (total volume = 25 μL). Incubate for 10 minutes at 4°C.
 - For mouse cells add 0.25 μL TruStain FcX™ PLUS (anti-mouse CD16/32) blocking reagent (BioLegend Cat. No. **156603**) to 5 x 10⁵ cells in 24.75 μL Cell Staining Buffer (total volume = 25 μL). Incubate for 10 minutes at 4°C.
- Transfer 25 μL of reconstituted cocktail to the tube containing 25 μL of FcR blocked cells. The final staining volume is 50 μL.
- Proceed with corresponding staining protocol:
 - For TotalSeq™-A panels, start at step 8 of the following protocol:
[go.biolegend.com/totalseq-a-v3-protocol](https://www.biolegend.com/totalseq-a-v3-protocol)
 - For TotalSeq™-B or TotalSeq™-C panels, start at step 8 of the following protocol:
[go.biolegend.com/totalseq-b-c-protocol](https://www.biolegend.com/totalseq-b-c-protocol)