

Tissue Dissociation Resources for use with the Milo Single-Cell Western System

The Milo Single-Cell Western system can analyze multiplexed protein expression patterns in dissociated tissues to identify and quantify cell subtypes contained in heterogeneous tissues. Generating high quality single-cell suspensions from tissues is an important precursor to a Single-Cell Western experiment in order to preserve cell states and relative abundance in downstream analysis. While we have not validated protocols for all tissue types, we hope these resources may be helpful when considering methods for dissociating your tissue of interest:

- **Demonstrated tissue dissociation protocol:** [Dissociation of Mouse Neural Tissue for Single-Cell Western Analysis](#)
- **Published methods:** Please refer to the [list of published Single-Cell Western papers](#) to see if there is a published protocol for your tissue of interest
- **Worthington Guide:** Please consult the [Worthington Tissue Dissociation Guide](#) which contains published tissue dissociation methods for a wide variety of tissue types
- **Commercial Options:** There are also several commercially available solutions for tissue dissociation methods including from Miltenyi Biotec, Sigma-Aldrich and Roche Diagnostics. Please consult their websites for further details.



Toll-free: (888) 607-9692
Tel: (408) 510-5500
info@proteinsimple.com
proteinsimple.com

For further assistance, please consult
support@proteinsimple.com

©2019 ProteinSimple. ProteinSimple, the ProteinSimple logo, Single-Cell Western and Milo are trademarks and/or registered trademarks of ProteinSimple.

PL9-0038, Rev A