

**ReadiUse™ TCEP removal solution**

Catalog number: 5540

Unit size: 1 ml

Component	Storage	Amount
-----------	---------	--------

**OVERVIEW****AT A GLANCE**

**Important** Warm up Readilink™ TCEP Removal Buffer at room temperature before use.

**SAMPLE EXPERIMENTAL PROTOCOL**

1. Estimate the TCEP amount to be removed.
2. Add the Readilink™ TCEP removal buffer to a solution that contains residual TCEP at molar ratio of 100 : 1 TCEP Removal Buffer/TCEP (to be removed).
3. Incubate the solution at room temperature from 30 to 60 minutes.
4. The solution is now ready to use for desired downstream assays.

**EXAMPLE DATA ANALYSIS AND FIGURES**

Example data analysis and images of this product can be found on the web at:  
<https://www.aatbio.com/products/readiuse-tcep-removal-buffer>

**DISCLAIMER**

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email [info@aatbio.com](mailto:info@aatbio.com) if you have any questions.