

**ReadiUse™ TCEP removal buffer**

Catalog number: 5540

Unit size: 1 ml

Component	Storage	Amount
ReadiUse™ TCEP removal buffer	Freeze (<-15 °C), Minimize light exposure	1 ml

**OVERVIEW**

TCEP is a common reducing reagent widely used to break disulfide bonds within and between proteins. Compared to the other two most common reducing agents, i.e., dithiothreitol (DTT) and 2-mercaptoethanol, TCEP has the advantages of being odorless, a more powerful reducing agent, an irreversible reducing agent, more hydrophilic, and more resistant to oxidation in air. TCEP is also used in the tissue homogenization process for RNA isolation. However, there are numerous reports that TCEP is not compatible with downstream assays that often have components reactive to TCEP. Our ReadiUse™ TCEP removing buffer can be effectively used for eliminating residual TCEP. It is a preformatted 2M solution that can be simply added to a solution that contains residual TCEP.

**DISCLAIMER**

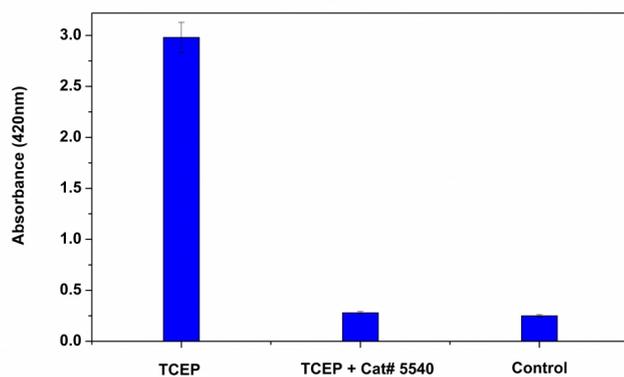
AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited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

**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**

**Figure 1.**

Reduction of Ellman's reagent by TCEP after it is treated with/without ReadiUse™ TCEP removal buffer (Cat #5540). TCEP was treated with/without ReadiUse™ TCEP removal buffer for 10min, and then the mixture was incubated with Ellman's reagent for 30min. Absorbance at 420nm was monitored using SpectraMax microplate reader.