

## 10X Protein Stabilization Solution

---

CATALOG NUMBER: PRB-5210

QUANTITY: 25 mL

### **Background**

Cell Biolabs' proprietary 10X Protein Stabilization Solution preserves long term activity of proteins during storage.

- Stabilizes and prevents aggregation of most enzymes and other proteins when compared to traditional methods such as glycerol
- Protects proteins from challenging environmental conditions such as high temperature fluctuation or freeze/thawing
- Solution contains low molecular weight components that can be removed downstream by dialysis or gel filtration

Cell Biolabs' proprietary 10X Protein Stabilization Solution is provided as an easily pipettable, buffered 10X concentrate which is added directly to the protein (do NOT pre-dilute). Subsequent to addition of the solution, proteins are stored at their normally recommended storage temperature from 4°C to -80°C. Although the level of stabilization is specific to each protein, the stabilizer maintains most protein activity compared to simply maintaining appropriate buffer and salt conditions. The 10X Protein Stabilization Solution is nontoxic, but solution components can be removed by dialysis or desalting before use in downstream assays, if desired.

### **Storage**

Store at -80°C. When ready to use, thaw and warm to 55°C, mix thoroughly to resuspend any precipitated material, and store on ice until use. Return the unused portion to -80°C.

### **Shelf Life**

One year from date of receipt under proper storage conditions.

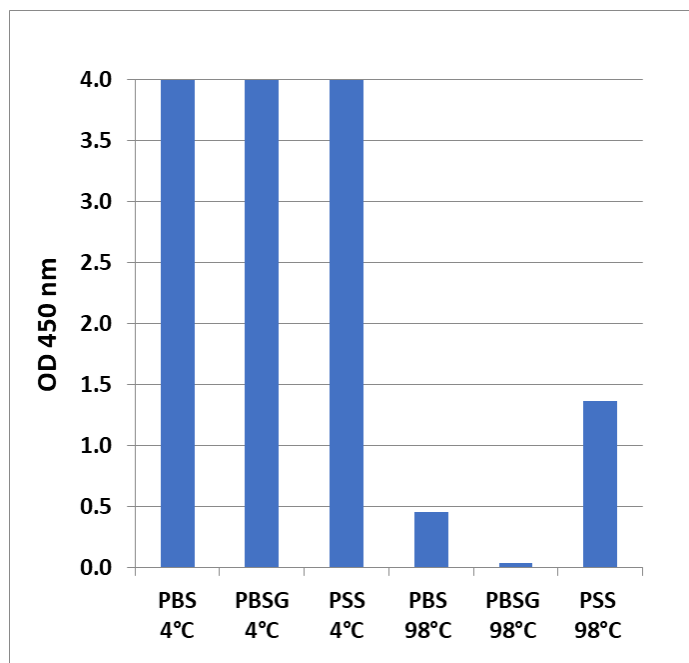
### **Protocol:**

1. Add 1/10 volume of 10X Protein Stabilization Solution to your protein of interest. Mix gently but thoroughly.
2. Store at the appropriate temperature for the protein of interest from 4°C to -80°C.

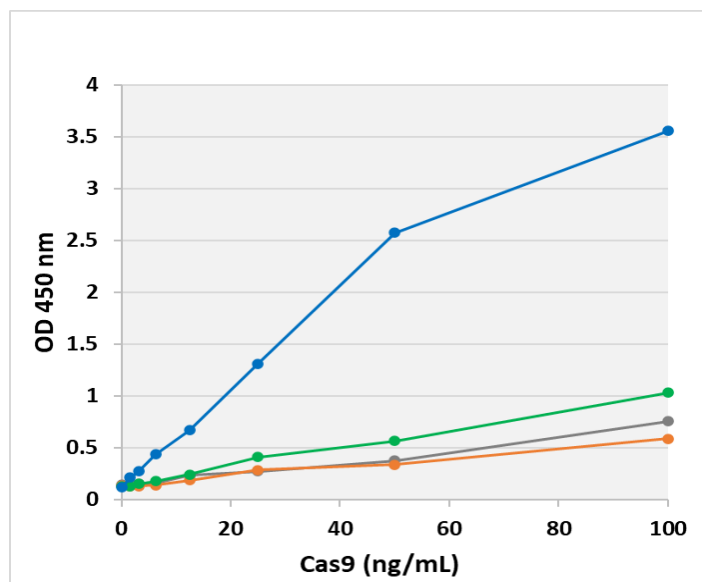
*Note: The 10X Protein Stabilization Solution does not change the storage properties of your protein of interest. Some proteins are inherently unstable at freezing temperatures and should be stored refrigerated.*

### Example of Results

The following figures demonstrate typical results. One should use the data below for reference only. This data should not be used to interpret actual results.



**Figure 1: Stabilization of Horseradish Peroxidase (HRP).** Purified HRP was incubated at 4°C or 98°C for 60 minutes and then tested for its activity by reaction with an ELISA Substrate Solution in the presence of phosphate buffered saline (PBS), PBS + 50% glycerol (PBSG), or PBS + Protein Stabilization Solution (PSS).



**Figure 2: Stabilization of Recombinant Cas9.** Purified recombinant Cas9 was incubated at 25°C for 6 days in 20 mM Tris pH 8, 300 mM NaCl, 10 mM EDTA containing 5% glycerol (grey line), 20%

glycerol (orange line), RIPA buffer (green line), or Protein Stabilization Solution (blue line), and then tested in the Cas9 ELISA kit (Cat No PRB-5079).

### **Warranty**

These products are warranted to perform as described in their labeling and in Cell Biolabs literature when used in accordance with their instructions. THERE ARE NO WARRANTIES THAT EXTEND BEYOND THIS EXPRESSED WARRANTY AND CELL BIOLABS DISCLAIMS ANY IMPLIED WARRANTY OF MERCHANTABILITY OR WARRANTY OF FITNESS FOR PARTICULAR PURPOSE. CELL BIOLABS' sole obligation and purchaser's exclusive remedy for breach of this warranty shall be, at the option of CELL BIOLABS, to repair or replace the products. In no event shall CELL BIOLABS be liable for any proximate, incidental or consequential damages in connection with the products.

***This product is for RESEARCH USE ONLY; not for use in diagnostic procedures.***

### **Contact Information**

Cell Biolabs, Inc.  
5628 Copley Drive  
San Diego, CA 92111  
Worldwide: +1 858-271-6500  
USA Toll-Free: 1-888-CBL-0505  
E-mail: [tech@cellbiolabs.com](mailto:tech@cellbiolabs.com)  
[www.cellbiolabs.com](http://www.cellbiolabs.com)

©2026: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.