

Stop Solution

Catalog Numbers SS03, SS04

Pub. No. MAN0017978 Rev. B (31)

Product description

Lot Number	See product label
Volume	25 mL (Cat. No. SS03), 1 L (Cat. No. SS04)
Product Description	ELISA Stop Solution is an acidic solution and should be handled with caution. ELISA Stop Solution is ready for use. Dilution before use is not recommended.
Applications	This solution is intended for use as a stop solution in enzyme assays which employs horseradish peroxidase and TMB substrate solution. It is suitable for use in microwell assays, such as ELISA.
Directions for Use	<p>The volume required for each assay should be decanted from the bottle into a clean dispensing trough. In microwell-based assays, the horseradish peroxidase in the wells is first incubated with TMB substrate solution. The volume of TMB substrate solution used for this application is usually 100 µL. The incubation time is usually 20–30 minutes. During incubation, the TMB substrate solution is converted from a colorless solution to a solution having a pronounced blue color as it reacts with the horseradish peroxidase. At the end of this incubation time, a volume of stop solution equal to the volume of TMB substrate solution is added to the wells to terminate the horseradish peroxidase reaction and the two solutions are allowed to mix in the wells.</p> <p>Addition of stop solution results in the rapid formation of a yellow end product. The optical density of this yellow end product is measured at 450 nm. The yellow end product is stable for sixty minutes after addition of stopping solution; therefore, the optical density should be measured within that time.</p>
Storage	Store at room temperature. It is highly recommended that the volume required for each assay be carefully decanted into a clean dispensing trough. The unused solution remaining in the trough should be discarded.
Expiration Date	See label

Limited product warranty

Life Technologies Corporation and its affiliates warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have questions, contact Life Technologies at www.thermofisher.com/support.



Bender MedSystems GmbH | Campus Vienna Biocenter 2 | 1030 Vienna, Austria

Revision history: Pub. No. MAN0017978 B (31)

Revision	Date	Description
B (31)	14 May 2024	Temperature factor changed under product description.
A (30)	10 July 2018	New document for ELISA Stop Solution - Product information sheet

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2018-2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.