



3050 Spruce Street
Saint Louis, Missouri 63103 USA
Telephone (800) 325-5832 (314) 771-5765
Fax (314) 286-7828
email: techserv@sial.com
sigma-aldrich.com

Product Information

Chemiluminescent Peroxidase Substrate for ELISA

Product Codes **CPS-2-60, CPS-2-120, and CPS-2-300**
Storage Temperature 2-8 °C

TECHNICAL BULLETIN

Product Description

Sigma's Chemiluminescent Peroxidase Substrate can be used for the highly sensitive detection of peroxidase conjugates in a variety of ELISA. This substrate is an enhanced luminol product with a stabilized peroxide buffer solution that provides a 100-fold improvement in signal to noise ratio over traditional luminol peroxidase formulations. The ultra high sensitivity combined with low background signal produces outstanding ELISA results.

Components

Chemiluminescent Peroxidase Substrate is available in 3 package sizes each containing Chemiluminescent Reagent (Product Code C 9986) and Chemiluminescent Reaction Buffer (Product Code C 0237).

Package Size	C 9986	C 0237
60 ml	20 ml	40 ml
120 ml	40 ml	80 ml
300 ml	100 ml	200 ml

Precautions and Disclaimer

This product is for laboratory research use only. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

Preparation Instructions

Prepare the Working Solution by mixing 1 part of the Chemiluminescent Reagent (C 9986) with 2 parts of the Chemiluminescent Reaction Buffer (C 0237). Mix well and protect from light.

Storage/Stability

The components are stable for a minimum of 18 months when stored at 2-8 °C in the original containers and protected from light. The Working Solution is stable for several hours at room temperature when protected from light.

Procedure

Sigma's Chemiluminescent Peroxidase Substrate is very sensitive and great care must be taken to optimize the individual assay components (antibodies, conjugates, etc).

For ELISA procedures, pipette 100 µl of the Working Solution into the appropriate wells, incubate for 5-10 minutes, and read using a luminometer (0.2-1 second integration time) or alternative method. Readings can be taken up to 60 minutes after addition of the Working Solution with minimal change in signal to noise ratios.

RBG,MAM 02/05-1

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.